

This version: US RUO E04A 070212 Version 2

of IFY is to be used only for reference purposes. Contents especially technical, methodical and regulations data, may not be amended.

IGFBP-3 LIA

Ligand-binding Immunoassay
for Quantitative Determination of

Functional Insulin-Like Growth-Factor Binding Protein-3

English

For Research Use Only.
Not for use in diagnostic procedures.

distributed in the US/Canada by:

EAGLE BIOSCIENCES, INC.

20A NW Blvd, Suite 112 Nashua, NH 03063

Phone: 617-419-2019 FAX: 617-419-1110

www.EagleBio.com • info@eaglebio.com



EAGLE
BIOSCIENCES



REF **E04A**

 **mediagnost**[®]

Gesellschaft für Forschung und Herstellung von Diagnostika GmbH



: Aspenhastr. 25 • D-72770 Reutlingen / Germany
Telefon: + 49 - (0) 7121 51484-0 • Fax: + 49 - (0) 7121 51484-10
E-Mail: contact@mediagnost.de • <http://www.mediagnost.de>

Symbols / Symbole

according to DIN EN 980 and EDMA recommendations Standard News 6 2001



Expiry date / Verfallsdatum / Date de péremption / Data di scadenza / Fecha de caducidad / Data de validade / Uiterste gebruiksdatum / Udløbsdato / Utgångsdatum / Termin Ważności / Lejárati idő / Doba expirácie / Doba expirace



Consider instructions for use / Bitte Gebrauchsanweisung beachten / Consultez la notice d'utilisation / Consultare le istruzioni per l'uso / Consulte las instrucciones de utilización / Consulte as instruções de utilização / Raadpleeg de gebruiksaanwijzing / Se brugsanvisningen / Läs anvisningarna före användning / Proszę przeczytać instrukcję wykonania / Vegyük figyelembe a használati utasításban foglaltakat / Postupujte podľa pokynov na použitie / Postupujte dle návodu k použití



Lot-Batch Number / Charge-Chargennummer / Lot-Code du lot / Lotto-Numero di lotto / Lote-Código de lote / Lote-Código do lote / Lot-Partijnummer / Lot-Batchkode / Lot-Satskod / Numer serii / Lot-Batch szám / Číslo šarže / Číslo šarže



Manufactured by / Hergestellt von / Fabriqué par / Prodotto da / Fabricado por / Fabricado por / Vervaardigd door / Fabrikation af / Tillverkad av / Wyprodukowane pr / Gyártotta / Vyrobené / Vyrobeno



Catalogue Number / Bestellnummer / Numéro de référence / Numero di riferimento / Número de referencia / Número de referência / Referentienummer / Referencenummer / Katalognummer / Numer katalogowy / Katalógusszám / Katalógové číslo / Katalogové číslo



Store at between / Lagerung bei zwischen / Conserver à entre / Conservare a tra / Conservar a temp. entre / Armazene a entre / Bewaar bij tussen / Opbevaars mellem / Förvaras vid / Przechowywać w / Tároljuk ő ő között / Skladujte v rozsahu / Skladujte v rozmezi



Contains sufficient for x tests / Inhalt ausreichend für x Tests / Contient suffisant pour x tests / Contenido suficiente per x test / Contiene suficiente para x pruebas / Contém suficiente para x testes / Bevat voldoende voor x bepalingen / Indeholder tilstrækkeligt til x prøver / Innehållet räcker till x analyser / Zawartość na x testów / Tartalma X teszt elvégzésére elegendő / Obsahuje materiál pre x testov / Obsahuje materiál pro x testů



Keep away from sunlight / Nicht dem Sonnenlicht aussetzen



Incubation time / Inkubationszeit



Incubate at / Inkubation bei



Shaking / schütteln



Mikrotiter plate/Mikrotiterplatte



Reconstitute in / Rekonstituieren in



Sample / Probe



LK Ligand Conjugate / Ligandkonjugat



EK Enzyme Conjugate / Enzymkonjugat



VP/PP Dilute in Buffer X / Verdünnen in Puffer X



A-E Standard X / Standard X



KS1/KS2 Control Serum / Kontrollserum



WP Washing Buffer Concentrate / Waschpufferkonzentrat



Washing Buffer / Waschpuffer



S Substrate / Substrat



SL Stop Solution / Stopplösung



Cover Plate with sealing tape / Platte abkleben



Measure plate within 30 min at 450 nm (Referencefilter ≥ 590 nm) / Ausmessung innerhalb von 30 min bei 450 nm (Referenzfilter ≥ 590 nm).

Symbols / Symbole	2
TECHNICAL PROPERTIES AND APPLICATIONS	4
Mediagnost IGFBP-3 LIA E04A	4
INTRODUCTION	4
Assay Characteristics	6
WARNINGS AND PRECAUTIONS	8
REAGENTS PROVIDED	9
MATERIALS REQUIRED BUT NOT PROVIDED	9
TECHNICAL NOTES	9
SPECIMEN COLLECTION, PREPARATION AND STORAGE	11
ASSAY PROCEDURE	11
QUALITY CONTROL	12
CALCULATION OF RESULTS	12
Establishing the Standard Curve	12
LIMITATIONS	13
EXEMPLARY VALUES	13
REFERENCES	15
SUMMARY – Functional IGFBP-3 LIA E04A	19
REF E04A International Test description	20

For in vitro use only!

For Research Use Only!

For professional use only!

CAUTION: Not for human or animal therapeutic or diagnostic use.

Read entire protocol before use!

TECHNICAL PROPERTIES AND APPLICATIONS

Mediagnost IGFBP-3 LIA E04A

- ◆ For research and professional use only!
- ◆ Quantitative determination of **functional bioactive (IGF-I binding) IGFBP-3: convenient “ELISA”, no sample pretreatment**
- ◆ solely **functional bioactive IGF-I-binding IGFBP-3** is being quantified, proteolytic fragments without binding affinity to the **natural ligand IGF-I** are not measured
- ◆ **Inter-Assay variation** of max. **6.8%**, **Intra-Assay variation** of max. **5.6%**
- ◆ reliable, fast and simple: performance like a conventional ELISA
- ◆ **Total IGFBP-3** integrates the GH secretory state over days, with stable serum levels due to absence of circadian variation a single measurement is highly informative for diagnosis of GH deficiency or GH excess: **correspondent measurement of functional IGFBP-3** reflects the respective degree of fragmentation of IGFBP-3 and thus allows further interpretation of the results
- ◆ **Direct Correlation** to the quantitative results of the immune reactive total IGFBP-3 from the respective Mediagnost Kits (E03A, IGF-R10, IGF-R11). In retrospect, or, in simultaneous determinations out of the same sample dilution!
- ◆ Small sample requirement, thus ideal for pediatric patients

INTRODUCTION

MEASURING IGFBP-3

All currently existing IGFBP-3 immunoassays use the binding of specific anti-IGFBP-3 antibodies for signal generation and thus IGFBP-3 quantification. The failure of differentiation between complete IGFBP-3 molecules and their respective fragments (derived physiologically due to the different proteases activities) is unavoidable in this system. Because one molecule IGFBP-3 can be cleaved in several fragments often false high quantitative values are measured. Based on this methodology it is not possible to differentiate between high IGFBP-3 levels in fact, or, a high degree of fragmentation. The incidental attempts to use monoclonal antibodies with a binding region represented only by the intact IGFBP-3 molecule are indirect, imprecise and insufficient. The activities of all effective proteases, which have different sites of action and therefore generate different kind of fragments, are disregarded.

The Mediagnost IGFBP-3 LIA however enables to determine the **real functional and effective bioactive IGFBP-3**, functional in terms of binding ability of the mainly interesting natural ligand, namely IGF-I !

The new test principle (patent pending DE19719001) uses anti-IGFBP-3 antibodies immobilized on the microtitre plate and biotinylated IGF-I as ligand. The IGFBP-3 of the sample is bound to the microtitre plate, and only concomitant bound, by IGFBP-3, biotinylated IGF-I gives the specific signal of the test. Therefore only functional IGFBP-3 is quantified. The patented test format ensures an easy and reliable performance, simply like a conventional ELISA kit. Advanced and time- as well as labor-intensive biochemical analysis (e.g., by size chromatography or Western Blots, etc.), which moreover only allows an estimation of concentrations, has become redundant.

TOTAL IGFBP-3

Insulin-like growth factors (IGF)-I and -II are bound to specific binding proteins (IGFBPs) in the circulation. To date, at least six binding proteins can be distinguished on the basis of their amino acid sequence. They are designated as IGFBP-1, IGFBP-2, ...IGFBP-6 (1). Lately the discovery of a new IGFBP-7 has been discussed (2). The predominating IGFBP in blood is IGFBP-3. In contrast to the other binding proteins, IGFBP-3 has the unique property to associate with an acid-labile non-binding subunit (ALS) after binding of either IGF-I or IGF-II (3-5). Most of the IGFBP-3 in plasma is present as the high molecular weight ternary complex, however, small amounts of free IGFBP-3 are also found (6, 7).

The development of specific immunoassays for IGFBP-3, those also recognizing the complete high molecular weight complex, provided new insights into its regulation (6-9). On the basis of these findings total serum IGFBP-3 has proved to be an additional useful test in the repertoire of diagnostic tools for evaluation of growth disorders (7, 8).

Patients with GH deficiency have subnormal total IGFBP-3 levels. In contrast, most of the small statured children with normal GH secretion have levels within the normal range (Figure 1).

The separation of these two groups is easy. A single measurement of the total IGFBP-3 concentration is sufficient for the diagnosis of GH deficiency with high accuracy (7, 18). In small statured children total IGFBP-3 levels rise to normal range within several days of GH administration and remain normal during continuous GH treatment (Figure 2). Therefore, total serum IGFBP-3 measurements are also suited for evaluating the potential of a patient to respond to GH and for GH therapy monitoring (19). In other patients of severe short stature, e.g. Ullrich-Turner syndrome or Silver-Russell syndrome, IGFBP-3 levels were found normal (8) reflecting normal GH secretion.

In normal tall children and adolescents without excessive GH secretion or in patients with Sotos syndrome, total IGFBP-3 levels are normal or slightly increased. In contrast, children with pituitary gigantism or adults with acromegaly have clearly elevated levels (Figure 3) (6, 15) that normalize on successful treatment. Therefore, total IGFBP-3 is also a useful parameter for the detection of excessive GH secretion and monitoring therapy efficacy. In precocious puberty, total IGFBP-3 levels are clearly increased by chronological age, whereas patients with premature thelarche have total IGFBP-3 levels in the upper normal range (15). Several factors besides GH influence total IGFBP-3 levels: age including sexual development, nutrition, hypothyroidism, diabetes mellitus, liver function and kidney function. Total IGFBP-3 levels are decreased by malnutrition, although less than IGF-I, in hypothyroidism, in diabetes mellitus and in hepatic failure (6-8), but are obviously increased in chronic renal failure (6, 10, 11). Measurement over 24 hours revealed constant circadian levels (12, 13). For clinical practice, the most important regulatory factor is GH. Single total IGFBP-3 measurements correlate significantly with the logarithm of the integrated spontaneous GH secretion (8, 14). In patients with GH deficiency, total IGFBP-3 levels are subnormal and increase gradually to within the normal range after several days of GH administration (7, 8). The slow response to GH and constant circadian levels during chronic daily application of GH (13) suggest that IGFBP-3 reflects the GH secretory state over days.

FRAGMENTED IGFBP-3

By proteolytical cleavage of the ternary complex of IGFBP-3, the physiological storage of IGF-I in circulation, IGF-I is released and subsequently able to bind to its cellular receptor.

IGFBP-3 can be cleaved by several proteases: Plasmin; PSA; MMPs; CathepsinD, Thrombin, gamma NGF.

Cleavage results not only in free IGF-I but also in different IGFBP-3 fragments. Dependent on the active protease, 22 cleavage sites are known, mostly located in the variable and the N-terminal region but some also in the C-terminal part. Fragments of about 30, 20 and 15 kDa can be generated by proteolysis. Their existence was proven by western ligand blotting. The

corresponding immunoblot demonstrated that not all fragments are able to bind IGF-I any more. So cleavage of IGFBP-3 results in at least partial loss of IGF-I affinity (20).

IGFBP-3 proteolysis can be found in a number of body fluids, like synovial fluid, amnion fluid, seminal fluid, interstitial fluid, peritoneal fluid, lymph and serum, of course. In all these body fluids different proteases can be activated resulting in a different fragmentation pattern of IGFBP-3 of different relevance for physiology (21).

Beside pregnancy, where nearly all serum IGFBP-3 is fragmented, several pathological conditions are known where the fragmentation level of IGFBP-3 is changed.

In certain previous studies for determining the degree of fragmentation of IGFBP-3 more sophisticated and laborious biochemical methods were involved. In these studies, in healthy individuals around 25 – 30% of serum total IGFBP-3 was found to be proteolyzed.

Assay Characteristics

It utilizes specific and high affinity antibodies for IGFBP-3, the antibodies are immobilized on the microtitre plate. The ligand, biotinylated IGF-I is pre-dispensed in excess into the needed wells. The sample is diluted outside within a special dilution buffer (Sample Buffer PP); all naturally bound IGFs are thereby released from their binding proteins. By adding an aliquot of such diluted sample with free IGFBP-3 to the ligand containing wells, the biotinylated IGF-I occupies all existing specific binding sites of the IGFBP-3. All IGFBP-3 molecules are bound afterwards to the microtitre plate, however by using a Streptavidin-Peroxydase-(POD)-Conjugate only the complexes of IGFBP-3/biotinylated IGF-I are involved in signal generation (in the closing POD-substrate reaction), thus, only functional IGFBP-3 is being quantified !

The standards of the LIA E04A are made of **native and functional human IGFBP-3** in concentrations of **0.4, 2, 6, 15 and 30 ng/ml**.

The **analytical sensitivity** of the LIA E04A has been determined with **0.18 ng/ml** (2 SD of zero standard in 16fold determination).

The Mediagnost IGFBP-3 LIA E04A is over a very wide range dilution true. The **Linearity of the dilution of sera** is excellent (s. Table 1).

Table 1: The linearity of the sample dilution

Dilution:	Sample 1 (recalculated, ng/ml)
1:100	2817
1:200	2962
1:400	3094
1:600	3055
1:800	3229
1:1000	2951
1:1200	2908
1:1600	2989
AV / SD / VC%	3001 / 125 / 4.2

AV = Average Value, **SD** = Standard deviation **VC** = Variation Coefficient%

The **Inter-** and **Intra-Assay** variation coefficients were found lower than **6.8% and 5.6%**. Exemplary determinations are shown in table 2 and table 3.

Table 2: Inter-Assay-Variation (n=8)

	Mean Value (ng/ml)	Standard Deviation (ng/ml)	VC (%)
Sample 1	1051	30	3.0
Sample 2	1891	129	6.8
Sample 3	2417	158	6.5

Table 3: Intra-Assay-Variation (n=16)

	Mean Value (ng/ml)	Standard Deviation (ng/ml)	VC (%)
Sample 1	1061	26	2,5
Sample 2	1571	85	5.4
Sample 3	2660	148	5.6

FRAGMENTED IGFBP-3

The results determined with the human IGFBP-3 LIA E04A are only for research use! Due to the new developed test system there are at present no concrete and specific clinical data available. Instructive might be the comparison of total IGFBP-3 values versus functional IGFBP-3 values in different subsets of samples.

Figure 4 shows results of comparative determinations of sera of healthy blood donors with the functional IGFBP-3 LIA E04A for functional IGFBP-3 and for total immunoreactive IGFBP-3 with Mediagnost kits, respectively. Functional IGFBP-3 values were found consistently lower on average compared with the corresponding total IGFBP-3.

Thereby a slight gender difference was obvious. Total IGFBP-3 serum concentrations of males were found lower than those of females. Functional IGFBP-3 concentrations of females and males however were found with nearly equal absolute values. Functional IGFBP-3 serum concentrations were found on average 25 % lower (Tab. 4; Fig 4).

Tab. 4: Mean total and functional IGFBP-3 values of sera of 103 female and 109 male healthy blood donors (IGFBP-3 concentrations in ng/ml). Average age was 42 years in each subset.

	Total IGFBP-3; min/max (ng/ml)	Functional IGFBP-3; min/max (ng/ml)	Functional IGFBP-3 (% of total)
Females	3568; 1752/5752	2506; 1102/4346	70,2
Males	2752; 1258/4726	2332; 1337/4304	84,8
overall	3160	2419	77,5

In a subset of pathological sera of donors suffering from different diseases substantially lower functional IGFBP-3 concentrations were obvious (Fig. 5). Further and more detailed studies however are necessary to reveal the underlying mechanisms and, from this to develop new diagnostic insights and processes.

WARNINGS AND PRECAUTIONS

The functional **IGFBP-3 LIA, E04A** is for in-vitro use only!. **For professional use only.**

Before starting the assay, read the instructions completely and carefully. Use the valid version of the package insert provided with the kit. Be sure that everything is understood. The Mediagnost GmbH is not liable for any loss or harm caused by non-observance of the instructions, as far as no law withstands.

Temperature WILL affect the absorbance readings of the assay. However, values for the patient samples will not be affected.

The shelf life of the components after opening is not affected, if used appropriately.

Do not use expired reagents.

Use separate pipette tips for each sample, control and reagent to avoid cross contamination. Use reservoirs only for single reagents. This especially applies to the substrate reservoirs. Using a reservoir for dispensing a substrate solution that had previously been used for the conjugate solution may turn solution colored. Do not pour reagents back into vials as reagent contamination may occur. Mix the contents of the microplate wells thoroughly to ensure good test results. Do not reuse microwells. Do not let wells dry during assay; add reagents immediately after completing the rinsing steps.

Caution: This kit contains material of human and/or animal origin.

Human Serum

Contained in following components: **Control Seren KS1 und KS2**

The source of human serum was tested by FDA recommended methods and found non-reactive for Hepatitis-B surface antigen (HBsAg), Hepatitis C virus (HCV), and Human Immunodeficiency Virus 1 and 2 (HIV) antibodies. No known test methods can offer total assurance of the absence of infectious agents; therefore all components and patient's specimens should be treated as potentially infectious.

Stop solution contains 0.2 M Sulphuric Acid (H₂SO₄)

R36/38 Irritating to eyes and skin

S26 In case of contact with eyes, rinse immediately with plenty of water and seek medical advice

S28.1 After contact with skin, wash immediately with plenty of water

S36/37 Wear suitable protective clothing and gloves

2-Methyl-4-Isothiazolin-3-one

contained in following components: **AK, VP, PP**

< 0.01% 2-Methyl-4-isothiazolin-3-one Solution

R34 Irritating to eyes and skin

R43 Sensibilisation through skin contact possible

S26 In case of contact with eyes, rinse immediately with plenty of water and seek medical advice

S36/37 Wear suitable protective clothing and gloves

S45 In case of accident or if you feel unwell seek medical advice

5-chloro-2-methyl 2H isothiazol-3-one and 2-methyl-2H-Isothiazol-3-one

contained in following components: **AK, VP, WP, PP**

< 0.01% (w/w) 5-chloro-2-methyl 2H isothiazol-3-one and 2-methyl-2H-Isothiazol-3-one Solution

R36/38 Irritating to eyes and skin

R43 Sensibilisation through skin contact possible

S26 In case of contact with eyes, rinse immediately with plenty of water and seek medical advice

S28.1 After contact with skin, wash immediately with plenty of water

TMB-Substrate (S) contains 3,3',5,5' Tetramethylbenzidine.

R20/21/R22 Harmful by inhalation, in contact with skin and if swallowed

R36/37/38 Irritating to eyes, respiratory system and skin

S26 In case of contact with eyes, rinse immediately with plenty of water and seek medical advice

S28.1 After contact with skin, wash immediately with plenty of water

S36/37 Wear suitable protective clothing and gloves

General first aid procedures:

Skin contact: Wash affected area thoroughly with water. Discard contaminated cloths and shoes.

Eye contact: In case of contact with eyes, rinse immediately with plenty of water at least 15 minutes.

In order to assure an effectual rinsing spread the eyelids.

Ingestion: If swallowed, wash out mouth thoroughly with water. Immediately see a physician.

Do not eat, drink or smoke in these areas.

Never pipette the materials with the mouth.

Spilled material must be wiped off immediately and should become disinfected. Clean contaminated areas and equipment with a suitable detergent.

REAGENTS PROVIDED

1)	MTP	Microtiter plate , ready for use, Microtiter plate with 96 wells, divided up in 12 strips à 8 wells (separately breakable). Coated with an antibody against human IGFBP-3.
2)	CAL	Standards A-E, 1 ml , lyophilised, contain native human functional IGFBP-3. Standard values are between 0.4 - 30 ng/ml (0.4, 2, 6, 15 and 30 ng/ml) functional IGFBP-3. Standards are to be reconstituted with 1 ml Sample Buffer PP each. Use 50 µl per well in the assay.
3)	DILU	Dilution Buffer VP, 30 ml , ready for use, please use for the dilution of the Ligand Conjugate LK and the Enzyme Conjugate EK
4)	DILU	Sample Buffer PP, 120 ml , ready for use, green colored, please use for reconstitution of Standards and Control and for dilution of the Samples and Control .
5)	Control	Control Seren KS1 and KS2, 250 µl , lyophilised, contain human Serum and should be reconstituted in 250 µl Sample Buffer PP. The functional IGFBP-3 target values and the respective ranges are given on the vial labels. The dilutions should be according to the dilution of the respected samples. Use 50 µl per well in the assay.
6)	Ligand	Ligand Conjugate LK, 100 µl , 101-fold Concentrate, contains biotinylated recombinant human IGF-I. Before use dilute 1:101 with VP. Use 50 µl per well in the assay.
7)	CONJ	Enzyme Conjugate EK, 140 µl , 101-fold Concentrate, contains HRP (Horseradish-Peroxidase)-labelled Streptavidin. Before use dilute 1:101 with VP. Use 100 µl per well in the assay.
8)	WASHBUF 20x	Washing Buffer (WP), 50 ml , 20 X concentrated solution. Washing buffer has to be diluted 1:20 with A.dest. or demineralised water before use (e.g. add the complete contents of the flask 50 ml into graduated flask and fill with A.dest to 1000 ml). Attention: After dilution, the Washing Buffer is only 4 weeks stable, please dilute only according to requirements.
9)	SUBST	Substrate (S), 12 ml , ready for use, horseradish-peroxidase-(HRP)-substrate, stabilised H ₂ O ₂ Tetramethylbencidine.
10)	H ₂ SO ₄	Stopping Solution (SL), 12 ml , ready for use, 0.2 M sulphuric acid, Caution acid!
11)		Sealing tape for covering of the microtiter plate, 2 x, adhesive.

MATERIALS REQUIRED BUT NOT PROVIDED

Precision pipettes and multichannel pipettes with disposable plastic tips
 Distilled or deionized water for dilution of the Washing Buffer (WP)
 Vortex-mixer
 Microtiter plate shaker (350 rpm)
 Microtiter plate washer (recommended)
 Micro plate reader ("ELISA-Reader") with filter for 450 and ≥590 nm
 Polyethylen PE/Polypropylen PP tubes for dilution of samples

TECHNICAL NOTES

The assay has to be conducted strictly according the test protocol herein.

Reagents with different lot numbers cannot be mixed. The microtiterplate and reagents are stable until the indicated expiry, if stored unopened and protected from sunlight at 2 – 8°C.

The shelf life of the components after opening is not affected, if used appropriately.

Bring all reagents to room temperature (20 - 25°C) before use. Possible precipitations in the buffers have to be resolved before usage by mixing and / or warming.

Incubation at room temperature means: 20-25°C

The incubation steps should be performed at mean rotation frequency of a particularly suitable microtitre plate shaker. We are recommending 350 rpm. Due to certain technical differences deviations may occur, in case the rotation frequency must become adjusted. Insufficient shaking may lead to inadequate mixing of the solutions and thereby to low optical densities, high variations and/or false values, excessive shaking may result in high optical densities and/or false values.

Proper washing is of basic importance for a secure, reliable and precise performance of the test. Incomplete washing is common and will adversely affect the test outcome. Possible consequences may be uncontrolled unspecific variations of measured optical densities, potentially leading to false results calculations of the examined samples. Effects like high background values or high variations may indicate washing problems.

All washing must be performed with the provided washing buffer diluted to usage concentration. Washing volume per washing cycle and well must be 300 µl at least.

The danger of handling with potentially infectious material must be taken into account.

When using an automatic microtitre plate washer, the respective instructions for use must be carefully followed. Device adjustments, e.g. for plate geometry and the provided washing parameters, must be performed. Dispensing and aspirating manifold must not scratch the inside well surface. Provisions must be made that the remaining fluid volume of every aspiration step is minimized. Following the last aspiration step of each washing cycle, this could be controlled, and possible remaining fluid could then be removed, by inverting the plate and repeatedly tapping it dry on non fuzzy absorbent tissue.

Manual washing is an adequate alternative option. Washing Buffer may be dispensed via a multistepper device, a multichannel pipette, or a squirt bottle. The fluid may be removed by dynamically swinging out the microtitre plate over a basin. If aspirating devices are used, care has to be taken that the inside well surface is not scratched. Subsequent to every single washing step, the remaining fluid should be removed by inverting the plate and repeatedly tapping it dry on non fuzzy absorbent tissue.

Standards and Control

The Standards **A – E** and **Control Sera KS1 and KS2** are reconstituted with the **Sample Buffer PP** provided in the Kit. It is recommended to keep the reconstituted reagents at room temperature for 15 minutes and then to mix them thoroughly but gently (no foam should result) with a Vortex mixer.

Reconstituted components (Standards (A – E) and Control Sera (KS1/KS2)) should be stored at 2-8°C for up to 1 week. If longer storage time is needed, store the components frozen at -20°C or below. Freezing extends the expiry at least 2 months. Avoid repeated freeze-thaw cycles. In case you plan to perform multiple independent IGFBP-3 determinations over a longer period with one kit, you should aliquot the components prior to freezing into suitable smaller volumes. This is strongly recommended.

Ligand and Enzyme Conjugate

Use the Dilution Buffer **VP** for the dilution of Ligand Conjugate **LK** and Enzyme Conjugate **EK** concentrates. The diluted solutions are only limited stable at 2-8°C.

Washing Buffer

The required volume of washing buffer is prepared by 1:20 dilution of the provided 20-fold concentrate with deionised water. The diluted Washing Buffer is stable for max. 4 weeks at 2-8°C.

Substrate Solution

The **Substrate Solution S**, stabilised H₂O₂-Tetramethylbencidine, is photosensitive – store and incubate in the dark.

Microtiterplate

Store the once unused microtiter strips and wells together with the desiccant in the tightly closed clip lock bag at 2-8°C use in the frame provided. The labelled expiry is not influenced in case of proper storage.

SPECIMEN COLLECTION, PREPARATION AND STORAGE

Serum samples as well as Heparin- and EDTA-Plasma samples are suited. Possible dilution of the sample by the anticoagulant must be considered. Citrate-Plasma samples are not recommended, values are reduced. A special external sample preparation prior to assay is not required.

Slight Hemolysis of the samples obviously doesn't disturb the determination. An external sample preparation prior to assay is not required (see below). Samples should be handled as recommended in general: collected and refrigerated as fast as possible. In case there will be a longer period (>24 hours) between the sample withdrawal and determination, store the undiluted samples frozen at -20°C or below in tightly closable plastic tubes. Avoid on principal repeated freeze-thaw cycles of serum/plasma (if required, please sub-aliquot) although functional IGFBP-3 levels were found to be unaffected by a few cycles, (3x) in our experiments. The high sensitivity of the assay allows the functional IGFBP-3 measurement in small sample volumes.

In most determinations (e.g., Serum- or Plasma samples and no extreme values expected) the dilution of **1:505 with Sample Buffer PP is suitable**, thus the respective covered **assay range is 0.2 to 15.15 mg/l**. Where required, depending on the expected IGFBP-3-values, the dilution with **Sample Buffer PP** can be higher or lower. The IGFBP-3 concentrations maybe completely different in body fluids of human origin other than serum or in cell culture supernatants.

Suggestion for dilution protocol:

Pipette **1 ml Sample Buffer PP** (green colored) in PE-/PP-Tubes (application of a multi-stepper is recommended in larger series), add **10 μl Serum- or Plasma** (dilution 1:101). Add 400 μl Sample Buffer **PP** in an other PE-/PP-tube and 100 μl of the thoroughly mixed first dilution (dilution 1:5). After mixing use **50 μl** of this 1:505 diluted solution within 1 hour **per determination** in the assay (pipetting control = blue coloring of the solution in the wells).

ASSAY PROCEDURE

NOTES: All determinations (Standards, Control Serum and samples) should be assayed in duplicate. For optimal results, accurate pipetting and adherence to the protocol are recommended.

When performing the assay, the Standards, Control Sera and the samples should be pipette as fast as possible (e.g. <15 minutes). To avoid distortions due to differences in incubation times, the **Enzyme Conjugate**, the **Substrate Solution S** as well as the **Stop Solution SL** should be added to the plate in the same order and in the same time interval each, respectively.

- 1) Please dilute the Ligand Conjugate LK 1:101 with **Dilution Buffer VP** (for use of the entire Microtitre Plate you may need 4.8 ml, thus you may dilute 60 μl of concentrated **LK** with 6 ml of buffer **VP**).
- 2) Add in every needed well **50 μl** of the already 1:101 diluted **Ligand Conjugate LK**
- 3) Add **50 μl Sample Buffer PP** in positions A1/2 (=blank wells!)
- 4) Pipette in positions B1/2 **50 μl** each **Standard A (0.4ng/ml)**,
pipette in positions C1/2 **50 μl** each **Standard B (2 ng/ml)**,
pipette in positions D1/2 **50 μl** each **Standard C (6 ng/ml)**,
pipette in positions E1/2 **50 μl** each **Standard D (15 ng/ml)**,
pipette in positions F1/2 **50 μl** each **Standard E (30 ng/ml)**.

To control the correct test accomplishment **50 μl** of the 1:505 (or in respective dilution rate of the sample) in Sample Buffer **PP** diluted **Control Sera KS1** and **KS2** can be pipetted in positions G1/2 and H1/H2.

Pipette **50 µl each** of the **diluted sample** (generally 1:505 diluted in Sample Buffer **PP**) in the rest of the wells, according to requirements. Please mix the dilutions immediately after sample addition and use within 60 minutes.

- 5) Cover the wells with the sealing tape and incubate the plate for **2 hours** at **room temperature** (shake at 350 rpm).
- 6) After incubation aspirate the contents of the wells and wash the wells **5 times** with **300 µl Washing Buffer WP**.
- 7) Following the last washing step, pipette **100 µl** of the **1:101 Enzyme Conjugate EK** in each well (for use of the entire Microtitre Plate you may need 9.6 ml, thus you may dilute 120 µl of concentrated **EK** with 12 ml of buffer **VP**).
- 8) Cover the wells with the sealing tape and incubate **1 hour** at **room temperature** (shake at 350 rpm).
- 9) After incubation wash the wells **5 times** with **Washing Buffer WP** as described in step 6
- 10) Pipette **100 µl of the TMB-Substrate solution S** in each well.
- 11) Incubate the plate for **30 Minutes in the dark** at **room temperature**.
- 12) After incubation pipette **100 µl Stop Solution SL** in each well.
- 13) Measure the absorbance **within 30 minutes at 450 nm (Reference filter ≥590 nm)**.

QUALITY CONTROL

Good laboratory practice requires that controls be run with each calibration curve. A statistically significant number of controls should be assayed to establish mean values and acceptable ranges to assure proper performance. The test results are only valid if the test has been performed following the instructions. Moreover the user must strictly adhere to the rules of GLP (Good Laboratory Practice) or other applicable federal, state or local standards/laws. All standards and kit controls must be found within the acceptable ranges as stated on the QC Certificate. If the criteria are not met, the run is not valid and should be repeated. Each laboratory should use known samples as further controls.

CALCULATION OF RESULTS

Establishing the Standard Curve

For the evaluation of the assay it is preconditioned that the absorbance values of the blank should be 0.25, these of standard E should be above 1.0.

Samples, which yield higher absorbance values than Standard E, are beyond the standard curve, for reliable determinations these samples should be tested anew with a higher dilution.

The standards provided contain the following concentrations of functional IGFBP-3:

Standard	A	B	C	D	E
ng/ml	0.4	2	6	15	30

- 1) Calculate the **mean absorbance** value for the blank from the duplicated determination (well A1/A2).
- 2) Subtract the mean absorbance of the blank from the mean absorbances of all other values.
- 3) Plot the standard concentrations on the x-axis versus the mean value of the absorbance of the standards on the y-axis.
- 4) Recommendation: Calculation of the standard curve should be done by using a computer program, because the curve is in general (without respective transformation) not ideally described by linear regression. **A higher-grade polynomial**, or **four**

parametric logistic (4-PL) curve fit or **non-linear regression** are usually suitable for the evaluation (as might be spline or point-to-point alignment in individual cases).

- 5) The functional IGFBP-3 concentration in ng/ml of the samples can be calculated by multiplication with the respective dilution factor, Division by 1000 converts the values in µg/ml or, equal mg/Litre (Example: a measured value was 6 ng/ml, Sample was 1:505 diluted: $6 \times 505 = 3030$ ng/ml, or 3,03 µg/ml or 3,03 mg/L, according the requested unit).

LIMITATIONS

IGFBP-3 levels are strongly dependent on GH secretion. However, a number of factors influence its plasma concentration and should be taken into account for appropriate interpretation. Plasma levels decrease during fasting (more than 1 day), in malnutrition, malabsorption, cachexia, impaired hepatic function, hypothyroidism, and diabetes mellitus. They may also be decreased in chronic inflammatory disease and malignancy. Levels are increased in states of impaired renal function and precocious puberty. In clinical situations with hyperprolactinemia or in patients with craniopharyngeoma, normal levels may be observed despite GH deficiency. In certain physiological (e.g. pregnancy) and pathological states, IGFBP-3 may be degraded to smaller molecular size compounds (16, 17) by specific proteases which affect IGFBP patterns.

EXEMPLARY VALUES

Exemplary values of fragmented IGFBP-3 in relation to total IGFBP-3 of healthy blood donors are shown in figure 1. It is demonstrated that about 25% of the IGFBP-3 is fragmented and thus could be detected by biotinylated IGF-I. In table 5 distribution of IGFBP-3 in healthy blood donors of different age and sex are shown.

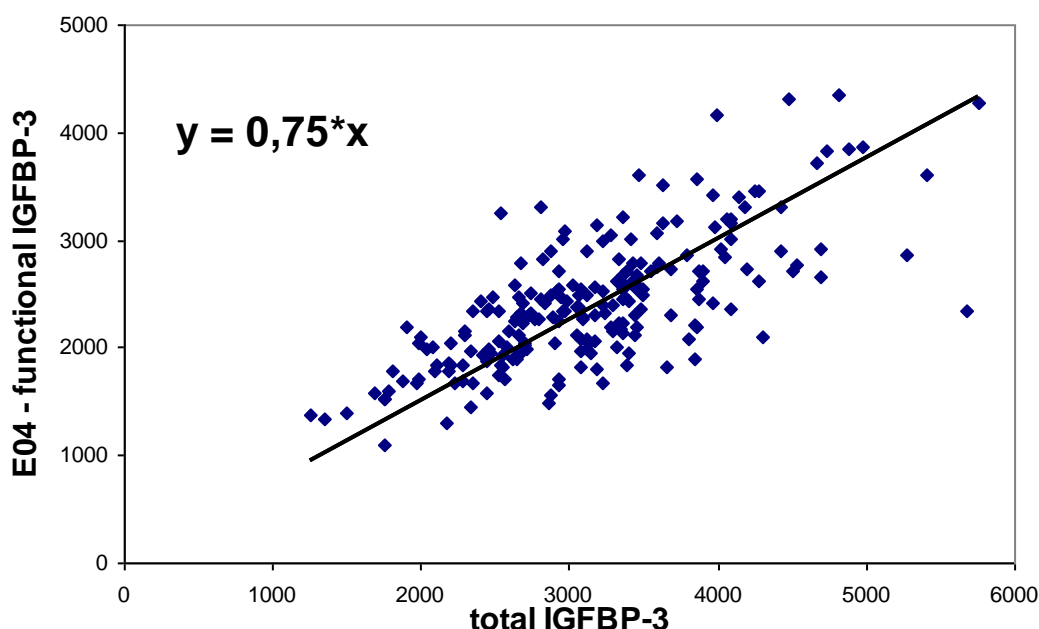


Figure 1: Plot of functional IGFBP-3 values against total IGFBP-3 values of sera from 212 healthy blood donors (IGFBP-3 concentrations in ng/ml).

Tab. 5: Serum levels of total immunoreactive IGFBP-3 in healthy subjects at various ages. Individuals between 7 and 17 years of age were classified according to gender, as the pubertal peak occurs almost 2 years earlier in girls than in boys.

Altersgruppe Age group	Percentiles / Perzentilen														
	0.1	1	5	10	20	30	40	50	60	70	80	90	95	99	
0-1 week	0.25	0.33	0.42	0.48	0.57	0.64	0.70	0.77	0.85	0.93	1.05	1.23	1.41	1.81	
1-4 weeks	0.49	0.62	0.77	0.86	0.99	1.10	1.19	1.29	1.40	1.52	1.68	1.93	2.16	2.68	
1-3 months	0.55	0.70	0.87	0.98	1.13	1.25	1.36	1.48	1.61	1.75	1.94	2.23	2.52	3.14	
3-6 months	0.64	0.80	0.98	1.10	1.25	1.38	1.49	1.61	1.74	1.88	2.07	2.37	2.65	3.24	
6-12 months	0.71	0.88	1.07	1.19	1.35	1.48	1.60	1.72	1.85	2.00	2.19	2.49	2.76	3.36	
1-3 years	1.02	1.21	1.41	1.53	1.69	1.82	1.94	2.05	2.17	2.31	2.48	2.74	2.98	3.47	
3-5 years	1.08	1.30	1.52	1.66	1.84	1.99	2.12	2.25	2.39	2.55	2.75	3.05	3.33	3.91	
5-7 years	1.19	1.42	1.66	1.81	2.01	2.16	2.30	2.44	2.59	2.76	2.97	3.29	3.59	4.2	
7-9 y.	boys	1.25	1.48	1.73	1.88	2.07	2.22	2.36	2.50	2.65	2.81	3.02	3.33	3.61	4.22
	girls	1.36	1.61	1.88	2.04	2.25	2.42	2.57	2.72	2.88	3.06	3.28	3.62	3.94	4.58
9-11 y.	boys	1.47	1.73	1.99	2.15	2.36	2.52	2.66	2.81	2.96	3.14	3.35	3.67	3.97	4.57
	girls	1.56	1.90	2.20	2.38	2.62	2.80	2.96	3.13	3.30	3.50	3.75	4.11	4.45	5.16
11-13 y.	boys	1.58	1.88	2.19	2.38	2.63	2.82	3.00	3.18	3.37	3.58	3.84	4.25	4.62	5.39
	girls	1.62	1.90	2.24	2.46	2.74	2.97	3.17	3.38	3.60	3.85	4.17	4.65	5.10	6.02
13-15 y.	boys	1.62	1.89	2.24	2.46	2.76	2.99	3.20	3.42	3.65	3.91	4.24	4.75	5.22	6.20
	girls	1.69	2.03	2.39	2.61	2.91	3.14	3.35	3.56	3.79	4.04	4.36	4.85	5.30	6.24
15-17 y.	boys	1.70	2.02	2.36	2.57	2.84	3.05	3.25	3.44	3.65	3.88	4.17	4.61	5.01	5.86
	girls	1.62	1.93	2.26	2.46	2.73	2.93	3.12	3.31	3.51	3.74	4.02	4.45	4.85	5.67
17-20 y.	1.58	1.90	2.24	2.45	2.72	2.94	3.13	3.33	3.54	3.78	4.07	4.53	4.95	5.83	
20-30 y.	1.55	1.86	2.20	2.41	2.68	2.90	3.09	3.29	3.50	3.74	4.04	4.50	4.92	5.80	
30-40 y.	1.44	1.75	2.08	2.29	2.56	2.78	2.98	3.18	3.39	3.64	3.95	4.42	4.86	5.78	
40-50 y.	1.38	1.68	2.01	2.21	2.48	2.69	2.88	3.08	3.29	3.53	3.83	4.29	4.72	5.63	
50-60 y.	1.34	1.64	1.96	2.16	2.42	2.63	2.83	3.02	3.23	3.46	3.76	4.22	4.65	5.55	
60-70 y.	1.28	1.58	1.90	2.10	2.37	2.58	2.78	2.98	3.19	3.44	3.75	4.23	4.67	5.62	
70-80 y	1.20	1.50	1.81	2.00	2.27	2.47	2.67	2.87	3.08	3.32	3.62	4.09	4.52	5.44	
> 80 y	1.13	1.43	1.73	1.92	2.19	2.39	2.59	2.79	3.00	3.23	3.54	4.00	4.44	5.36	

Serum levels are given as mg/L
 Die Serumkonzentrationen sind in mg/L angegeben
 week = Woche; months = Monate
 y. = years; =Jahre

Determined with IGFBP-3 RIA (Blum et al. 1990)
 Mit IGFBP-3-RIA gemessen (Blum et al. 1990)
 The values above 70 years are extrapolated.
 Die Werte für über 70-Jährige sind extrapoliert.

REFERENCES

- 1) Ballard J, Baxter R, Binoux M, Clemmons D, Drop S, Hall K, Hintz R, Rechler M, Rutanen E, Schwander J (1989) On the nomenclature of the IGF binding proteins. *Acta Endocrinol (Copenh)* 121:751-752
- 2) Wilson EM, Oh Y, Rosenfeld RG (1997) Generation and characterization of an IGFBP-7 antibody: Identification of 31 kD IGFBP-7 in human biological fluids and Hs578T human breast cancer conditioned media. *J Clin Endocrinol Metab* Vol 82, 4:1301-1303
- 3) Baxter RC (1988) Characterization of the acid-labile subunit of the growth hormone-dependent insulin-like growth factor binding protein complex. *J Clin Endocrinol Metab* 67:265-272
- 4) Baxter RC, Martin JL (1989) Structure of the Mr 140,000 growth hormone dependent insulin-like growth factor binding protein complex: determination by reconstitution and affinity-labeling. *Proc Natl Acad Sci USA* 86:6898-6902
- 5) Holman SR, Baxter RC (1996) Insulin-like growth factor-binding protein-3: factors affecting binary and ternary complex formation. *Growth Regulation* 6: 42-47.
- 6) Baxter RC, Martin J (1986): Radioimmunoassay of growth hormone-dependent insulin-like growth factor binding protein in human plasma. *J Clin Invest* 78:1504-1512
- 7) Blum WF, Ranke MB, Kietzmann K, Gauggel E, Zeissel HJ, Bierich JR (1990) A specific radioimmunoassay for the growth hormone (GH)-dependent somatomedin-binding protein: its use for diagnosis of GH deficiency. *J Clin Endocrinol Metab* 70:1292-1298
- 8) Blum WF, Ranke MB (1990) Use of insulin-like growth factor binding protein 3 for the evaluation of growth disorders. *Horm Res* 34 (Suppl):31-37
- 9) Blum WF (1993) Insulin-like growth factor-binding protein 3: Entwicklung eines Radioimmunoassays und Untersuchungen zur klinischen Bedeutung. Habilitationsschrift, Tübingen.
- 10) PDK, Hintz RL, Sperry JB, Baxter RC, Pow Fig. 7 1989) IGF-binding proteins in growth-retarded men with chronic renal failure. *Pediatr Res* 26:30
- 11) Blum WF, Ranke MB, Kietzmann K, Tönshoff B, Menis U (1989) Excess of-IGF-binding proteins in chronic renal failure: evidence for relative GH resistance and inhibition of somatomedin activity. In: Drop SLS, Hintz RL (eds) *Insulin-like Growth Factor Binding Proteins*. Excerpta Medica, Amsterdam, pp 93-101
- 12) Baxter RC, Cowell CT (1987) Diurnal rhythm of growth hormone-independent binding protein for insulin-like growth factors in human plasma. *J Clin Endocrinol Metab* 65:432-440
- 13) Jorgensen JOL, Blum WF, Moller N, Ranke MB, Christiansen JS (1990) Circadian patterns of serum insulin-like growth factor (IGF)-II and IGF-binding protein 3 in growth hormone deficient patients and age- and sex-matched normal subjects. *Acta Endocrinol (Copenh.)* 123:257-262
- 14) Blum WF, Albertsson-Wikland K, Rosberg S, Jorgensen JOL, Ranke MB (1990) Insulin-like growth factor binding protein 3 (IGFBP-3) reflects spontaneous growth hormone (GH) secretion. *Horm Res* 33 (Suppl 3): 3(abstract)
- 15) Blum WF, Ranke MB (1990) Insulin-like growth factor-binding proteins (IGFBPs) with special reference to IGFBP-3. *Acta Paediatr Scand (Suppl)* 367:55-62
- 16) Giudice LC, Farrell EM, Pham H, Lamson G, Rosenfeld RG (1990) Insulin-like growth factor binding proteins in maternal serum throughout gestation and in the puerperium: effects of a pregnancy-associated serum protease activity. *J Clin Endocrinol Metab* 71:806-816
- 17) Hossenlopp P, Segovia B, Lassarre C, Roghani M, Bredon M, Binoux M (1990) Evidence of enzymatic degradation of insulin-like growth factor-binding proteins in the 150k complex during pregnancy. *J Clin Endocrinol Metab* 71:797-805
- 18) Ranke MB, Schweizer R, Elmlinger MW, Weber K, Binder G, Schwarze CP, Wollmann HA (2000) Significance of Basal IGF-I, IGFBP-3 and IGFBP-2 Measurements in the diagnostics of short stature in children. *Horm Res* 2000;54:60-68
- 19) Ranke MB, Schweizer R, Elmlinger MW, Weber K, Binder G, Schwarze CP, Wollmann HA (2001) Relevance of IGF-I, IGFBP-3, and IGFBP-2 Measurements during GH treatment of GH-deficient and non-GH-deficient children and adolescents. *Horm Res* 2001; 55:115-124
- 20) Binoux M, L.C., Mohseni-Zadeh S. Biological Actions of Proteolytic Fragments of the IGF Binding Proteins. in *The IGF System* (ed. Rosenfeld R, R.C.) (Humana Press Inc, Totowa NJ, 1999).
- 21) Maile LA, C.A., Holly JMP. IGF Binding Protein Prtealysis in Various Clinical States. in *The IGF System* (ed. Rosenfeld R, R.C.) 633-649 (Humana Press Inc, Totowa, NJ, 1999).
- 22) Fielder PJ, G.-A.J., Rosenbloom AL, Carlsson L, Hintz RL, Rosenfeld RG. Expression of serum insulin-like growth factors, insulin-like growth factor binding proteins, and the growth hormone binding protein in heterozygote relatives of Ecuadorian growth hormone receptor deficient patients. *J Clin Endocrinol Metab* 74, 743-750 (1992).
- 23) Cotterill AM, H.J., Taylor AM, Davies SC, Coulson VJ, Preece MA, Wass JAH, Savage MO. The insulin-like growth factor (IGF-)binding proteins and IGF bioactivity in Laron-type dwarfism. *J Clin Endocrinol Metab* 74, 56-63 (1992).
- 24) Davies DC, W.J., Ross RJM, Cotterill AM, Buchanan CR, Coulson VJ, Holl JMP. The induction of a specific protease for insulin-like growth factor binding protein-3 in the circulation during severe illness. *J Endocrinol* 130, 469-473 (1991).

- 25) Cwyfan Hughes SC, C.A., Molloy AR, Cassel TMB, Braude N, Hinds CH, Wass JAH. The induction of specific proteases for insulin-like growth factor binding proteins following major heart surgery. *J Endocrinol* 135, 135-145 (1992).
- 26) Timmins AC, C.A., Hughes SCC, HOLly JMP, Ross RJM, Blum W, Hinds CJ. Critical illness is associated with low circulating concentrations of insulin-like growth factor I and growth factor II, alterations in insulin-like growth factor binding -proteins and induction of an insulin -like growth factor binding-protein-3 protease-. *Crit Care Med* 24, 1460-1466 (1996).
- 27) Bang, P., Brismar K, Rosenfeld RG, Hall K. Fasting affects serum insulin-like growth factors (IGFs) and IGF binding proteins differently in patients with non-insulin dependent diabetes versus healthy non-obese and obese subjects. *J Clin Endocrinol Metab* 78, 960-967 (1994).
- 28) Lang CH, F., Frost RA, Gelato MC, Sakurai Y, Herndon DN, Wolfe RR. Regulation of the insulin-like growth factor system by insulin burn patients. *J Clin Endocrinol Metab* 81, 2474-2480 (1996).
- 29) Frost VJ, H.S., Lonning PE, van der Stappen J, HOLly JMP. Effects of treatment with megestrol acetate, aminoglutethimide or formestane on insulin-like growth factor (IGF) I and II and IGF-binding proteins (IGFBPs) and IGFBP-3 protease status in patients with advanced breast cancer. *J Clin Endocrinol Metab* 81, 2216-2221 (1996).
- 30) Renehan AG, K.K., Diamandi A, O'Dwyer ST, Shalet SM. Quantification of serum IGFBP-3 proteolysis in patients with colorectal cancer. *Growth Horm IGF Res* 13, 217 (2003).
- 31) Muller HL, O.Y., Gargosky SE, Lehrnbecker T, Hintz RL, Rosenfeld RG. Concentrations of insulin-like growth factor (IGF) binding protein-3 (IGFBP-3), IGF and IGFBP-3 protease activity in cerebrospinal fluid of children with leukemia, central nervous system tumor or meningitis. *J Clin Endocrinol Metab* 77, 113-1119 (1993).
- 32) Fernihough JK, B.M., Cwyfan-Hughes S, Holly JMP. Local disruption of the insulin-like growth factor system in the arthritic joint. *Arthritis Rheum* 39, 1556-1565 (Local disruption of the insulin-like growth factor system in the arthritic joint.).
- 33) Lassarre C, M. Binoux. Measurement of intact Insulin-Like Growth Factor-Binding Protein-3 in human plasma using a ligand immunofunctional assay. *J Clin Endocrinol Metab* 86, 1260-1266 (2001)
- 34) Lassarre C, F. Duron, M. Binoux. Use of the Ligand Immunofunctional Assay for human Insulin-like growth factor (IGF) binding protein-3 (IGFBP-3) to analyze IGFBP-3 proteolysis and IGF-I bioavailability in Healthy adults, GH-deficient and acromegalic patients, and diabetics. *J Clin Endocrinol Metab* 86, 1942-1952 (2001)
- 35) Langkamp M. Weber K., Kirschner M., Pridzun L., Ranke M.B. Validation of Functional insulin-Like Growth Factor Binding Protein-3 measurement by a Ligand Immunoassay. *Clin Lab.* 11+12/2010, 56; 535-542

SUMMARY – Functional IGFBP-3 LIA E04A

Reconstitution/ Dilution of Reagents:		
Standards A-E	Reconstitution in Sample Buffer PP (green)	1 ml each
Control Sera KS1 & KS2	Reconstitution in Sample Buffer PP (green)	250 µl
Ligand Conjugate LK	Dilute LK and EK 1:101 each in Dilution Buffer VP (e.g. 60 µl LK plus 6 ml VP each)	1:101 each
Enzyme Conjugate EK	(e.g. 120 µl EK plus 12 ml VP each)	
Washing Buffer WP	dilute in A. dest. (e.g. add the complete contents of the flask 50 ml into a graduated flask and fill with A.dest. to 1000 ml)	1:20
Sample Dilution + Control Sera KS1 & KS2: 1:505 in Sample Buffer PP (green colored), mix directly and use within max. 60 min.		
Use 50 µl per determination (pipetting control= blue coloration)		
Before assay procedure bring all reagents to room temperature		

Proposal of Assay Procedure for Double Determination:

Pipette	Reagents	Well Positions
50 µl	diluted Ligand Conjugate LK	Pipette in <u>all</u> required number of wells
50 µl	Sample Buffer PP as Blank	A1 and A2
50 µl	Standard A (0.4 ng/ml)	B1 and B2
50 µl	Standard B (2 ng/ml)	C1 and C2
50 µl	Standard C (6 ng/ml)	D1 and D2
50 µl	Standard D (15 ng/ml)	E1 and E2
50 µl	Standard E (30 ng/ml)	F1 and F2
50 µl	diluted Control Serum KS1	G1 and G2
50 µl	diluted Control Serum KS2	H1 and H2
50 µl	diluted Samples	Pipette samples in the rest of the wells according to requirements
Cover the wells with the sealing tape		

Incubation: 2 h at RT, 350 rpm

5x 300 µl	Aspirate the contents of the wells and wash 5x with 300 µl each WP/ well	each well
100 µl	diluted Enzyme Conjugate EK	each well

Incubation: 1 h at RT, 350 rpm

5x 300 µl	Aspirate the contents of the wells and wash 5x with 300 µl each WP/ well	each well
100 µl	Substrate Solution S	each well


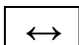
Incubation: 30 min in the dark at RT

100 µl	Stop Solution SL	each well
Measure the absorbance within 30 min at 450 nm (≥590 nm Reference)		


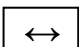
CAL A-E	STD A -E	Rec in 1ml PP	
Control	KS1 & KS2	Rec in 250 µl PP	1:505 DILU PP
LIGAND	LK		1:101 DILU VP
CONJ	EK		1:101 DILU VP
WASHBUF 20x	WP		1:20 DILU A. dest.

SPE	1:505 DILU PP
°C 20-25 °C	

50 µl	LK (1:101 DILU VP)	A1 - End
50 µl	PP	A1/2
50 µl	CAL A (0.4 ng/ml)	B1/2
50 µl	CAL B (2 ng/ml)	C1/2
50 µl	CAL C (6 ng/ml)	D1/2
50 µl	CAL D (15 ng/ml)	E1/2
50 µl	CAL E (30 ng/ml)	F1/2
50 µl	CONTROL KS1 1:505 DILU PP	G1/2
50 µl	CONTROL KS2 1:505 DILU PP	H1/2
50 µl	SPE 1:505 DILU PP	
TAPE		

 2 h **°C** 20-25  350 rpm

5x300 µl	5x WASHBUF WP
100 µl	EK (1:101 DILU VP)
TAPE	

 1 h **°C** 20-25  350 rpm

5x300 µl	5x WASHBUF WP
100 µl	SUBST TMB S

 0.5 h **°C** 20-25 

100 µl	H₂SO₄ SL
MEASURE	