

Mouse/Rat IGFBP-3 ELISA

Enzyme Immunoassay for Quantitative Determination of
Mouse- and Rat-
Insulin-like Growth Factor Binding Protein-3 (IGFBP-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

E031



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>

| | |
|---|--|
|  | Expiry date/ Verfallsdatum/ Date de péremption/ Data di scadenza/ Fecha de caducidad/ Data de validade/ Uiterste gebruiksdatum/ Udløbsdato/ Bäst före-datum/ Termin ważności/ Lejárati idő/ Țas expirăcie/ Doba expirace/ Срок на годност/ Aegumiskuupäev/ Ημερομηνία λήξης/ Data de expirare/ Rok uporabe/ Viimeinen käyttöpäivä |
|  | Consider instructions for use/ Bitte Gebrauchsanweisung beachten/ Consultez la notice d'utilisation/ Consultare le istruzioni per l'uso/ Consulte las instrucciones de uso/ Respeitar as instruções de utilização/ A.u.b de gebruiksaanwijzing volgen/ Se brugsanvisningen/ Läs anvisningarna före användning/ Proszę przeczytać instrukcję obsługi/ Vegye figyelembe a használati utasításban foglaltakat/ Postupujte podľa pokynov na použitie/ Dodržujte návod k použití/ Моля, спазвайте инструкцията за употреба/ Palun järgige kasutusjuhendit/ Λάβετε υπόψη σας τις οδηγίες χρήσης/ Vă rugăm să respectați instrucțiunile de utilizare/ Upošteevajte navodila za uporabo/ Lue käyttöohje huolellisesti! |
| g | 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/ Partisatskod/ Numer serii/ Tétel-sarzs szám/ Číslo šarže/ Číslo šarže/ Партиден номер/Partii number/ Παρτίδα-αριθμός παρτίδας/ Lot-număr lot/ Številka serije/ Erä |
| M | Manufactured by/ Hergestellt von/ Fabriqué par/ Prodotto da/ Fabricado por/ Fabricado por/ Vervaardigd door/ Fabrikation af/ Tillverkad av/ Wyprodukowane przez/ Gyártotta/ Vyrobene/ Vyrobeno v/ Производител/ Tootja/ Κατασκευάζεται από/ Produs de/ Proizvajalec/ Valmistaja |
| h | Catalogue Number/ Bestellnummer/ Numéro de référence/ Numero di riferimento/ Número de referencia/ Número de Referencia/ Referentienummer/ Referencenummer/ Beställningsnummer/ Numer katalogowy/ Rendelési szám/ Katalógovné číslo/ Objednací číslo/ Каталоген номер/ Tellimisnumber/ Αρ. παραγγελίας/Număr comandă/ Številka naročila/ Viite tai tilausnumero |
|  | Store at between/ Lagerung bei zwischen/ Conserver à entre/ Conservare a tra/ Conservar a temp. entre/ Armazena entre/ Bewaar bij tussen/ Opbevares mellem/ Förvaras vid/ Przechowywać w/ Tárolási tartomány/ Skladujte v rozsahu / Skladujte v rozmezí/ Температурно ограничение/ Säilitada temperatuuridel/ Φύλαξη σε θερμοκρασία/ Depozitare între/ Skladiščenje med/ Säilytys x-y Celsiusasteen lämpötilassa |
| X | Contains sufficient for x tests/ Inhalt ausreichend für x Tests/ Contient suffisant pour x tests/ Contenuto sufficiente per x test/ Contenido suficiente para x pruebas/ Conteúdo 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/ Obsah dostačuje pro x testů/ Съдържа достатъчно за x тестове/ Sisust jätkub x katse jaoks/ Το περιεχόμενο επαρκεί για x δοκιμές/ Conținut suficient pentru x teste/ Vsebina zadostuje za x preizkusov/ Sisältö riittää x testille |
|  | Keep away from sunlight/ Nicht dem Sonnenlicht aussetzen/ Conserver à l'abri de la lumière/ Conservare al riparo della luce solare/ No exponer a la luz solar/ Proteger da luz solar/ Niet aan zonlicht blootstellen/ Må ikke udsættes for sollys/ Utsatt inte för solljus/ Nie wystawiać na słońce/ Napfénytől távol tartandó/ Nevystavovat slnečnému svetlu/ Nevystavovat slunečnému světlu/ Да се предпазва от слънчева светлина/ Kaitsta otsese päikesekiirguse eest/ Κρατήστε το μακριά από την ηλιακή ακτινοβολία/ Τηνετή departe de lumina soarelui/ Ne izpostavljajte sončni svetlobi/ suojaa auringonvalolta |
|  | Incubation time/ Inkubationszeit/ Temps d'incubation/ Tempo d'incubazione/ Tiempo de incubación/ Tempo de incubação/ incubatietijd/ Inkubationstid/ inkubationstid/ Czas inkubacji/ Inkubációs idő/ Inkubačná lehota/ Inkubační doba/ Инкубационен период/ Inkubatsiooniaeg/ Χρόνος επώασης/ Timp de incubare/ Inkubacijska doba/ inkubaatioaika |
|  | Incubate at/ Inkubation bei/ Incuber à/ Incubare a/ incubar a/ Incubar a/ incubatietemperatuur/ Inkubation ved/ inkubation vid/ Inkubacja przy/ Inkubáció hőmérséklete/ Inkubácia pri/ Inkubace při/ Инкубира се при/ Inkubatsioon temperatuuril/ Επώαση στους/ Incubare la/ Inkubacija pri/ Inkubaatiolämpötila |
|  | Mix tubes with a Vortex mixer/ Mix Röhrchen mit Vortex Mixer/ Mélanger à l'aide d'un vortex/ Miscelare la provetta con agitatore Vortex/ Tubos de mezcla con mezclador de vortex/ Misturar os tubos com um agitador Vortex/ buisjes mengen met een Vortex/ Blanderør med Vortex-mixer/ Blanda rören med en vortexblandare/ Miksowanie rurek w mikserze Vortex/ Csővecskék keverése örvénykeverővel/ Premiešat pomocou prístroja Vortex/ Promíchat pomocí přístroje Vortex/ Разбъркване на епруветките с миксер Vortex/ Segada torukesi Vortexi mikseriga/ Αναμίξτε τους σωληνίσκους με αναδευτήρα Vortex/ Amestecați erubetele cu ajutorul unui agitator vortex/ Mešanje cevčic z mešalnikom Vortex/ Sekoita putket Vortex sekoittajalla |
| MTP | Mikrotiterplate/ Mikrotiterplatte/ plaque de microtitrage/ Piastra di microtitolazione/ Placa de microtitulación/ Placa de Microtitulação/ Microtiterplaat/ Mikrotiterplade/ mikrotiterplatta/ microtiterplaat/ Plytka microtiter/ Mikrotiter lap/ Mikrotitračná podložka/ Mikrotitrační podložka/ Микротитърна плака/ Mikrotiterplaat/ Τρυβλίο μικροπιλοδότησης/ Microplacă/ Mikrotitrská plošča/ Mikrotitrauslevy |
| Rec in | Reconstitute in/ Rekonstituieren in/ Reconstituer dans/ Ricostituire nel/ Reconstituir en/ Reconstituir em/ Reconstituieren in/ Rekonstituér i/ Rekonstituera/ Rekonstytiować w/ Helyreállítás/ Znovu pripravít za/ Znovu pripravít za/ Разтваряне в/ Moodustada uuesti/ Ανασυστήστε σε/ Reconstituire în/ Predelava v/ Rekonstituoi |
| SPE | Sample/ Probe/ Echantillon/ Campione/ Muestra/ Amostra/ Monster/ Prøve/ prov/ Próbka/ Minta/ Vzorka/ Vzorek/ Проба/ Proov/ Δείγμα/ Probă/ Vzorec/ Näyte |
| Ab | Antibody Conjugate/ Antikörperkonjugat/ Anticorps conjugué/ Coniugato di anticorpo/ Conjugado de anticuerpos/ Conjugado anticorpo/ Antilichaamconjugaat/ Antistoffer-konjugat/ Antikroppskonjugat/ Koniugat antyciał/ Antitest páros/ Protílátkový konjugát/ Protílátkový konjugát/ Антитяло конюгат/ Antikehad konjugaat/ Σύμπλοκο αντισώματος/ Compuși din anticorpi/ Antitelesa konjugat/ Vasta-aine konjugaati |

| | |
|---|--|
| CONJ | Enzyme Conjugate/ Enzymkonjugat/ Conjugué enzymatique/ Coniugato di enzima/ Conjugado de enzimas/ Conjugado Enzima/ Enzymkonjugaat/ Enzym-konjugat/ Enzymkonjugat/ Koniugat enzymów/ Enzim páros/ Enzymatický konjugát/ Enzymatický konjugát/ ензим конюгат/ Ensüümi konjugaat/ Σύμπλοκο –ενζύμου/ Compuși din enzime/ Encima konjugat/ Enzymi-konjugaatti |
| BUF | Buffer/ Puffer/ Tampon/ Tampone/ Tampón/ Tampão/ Buffer/ Buffer/ Buffer/ Bufor/ Puffer/ Pufer/ Pufri/ Буфер/ Puhver/ Ρυθμιστικό διάλυμα/ Tampon/ Puffer/ Puskuri |
| DILU BUF X | Dilute in Buffer X/ Verdünnen in Puffer X/ Diluer dans le tampon X/ Diluire nel tampone X/ Diluir en tampón X/ Diluir no Tampão X/ Verdunnen in buffer X/ Fortyndes i buffer X/ Späd i buffert X/ Rozcieńczenie w buforze X/ Hígítás X pufferben/ Riedit' v pufri X/ Ředit v pufru X/ Разреждане в буфер X/ Lahjendada puhvris X/ Αραιώστε σε ρυθμιστικό διάλυμα X/ Diluați în tamponul X/ Razredčiti v pufru X/ Laimennetaan x puskuriin |
| STD | Standard X/ Standard X/ Etalon X/ Standard X/ Estándar X/ Standard X/ Standaard X/ Standard X/ standard X/ Standard X/ Standard X/ Štandard X/ Standard X/ Стандарт X/ Standard X/ Πρότυπο X/ Standard X/ Standardni X/ Standardi X |
| Control | Control Serum X/ Kontrollserum X/ Contôle sérique X/ Siero di controllo X/ Suero de control X/ Soro de Controlo X/ controleserum X/ Kontrolserum X/ Kontrollserum X/ Serum kontrolne X/ Ellenőrző szérum X/ Kontrolné serum X/ Kontrolní serum X/ Контролен серум X/ Kontrollseerum X/ Ορός ελέγχου X/ Ser de control X/ Kontrolni serum X/ Kontrolli seerumi X |
| WASHBUF 20x | Washing Buffer Concentrate/ Waschpufferkonzentrat/ Tampon de lavage conc./ Tampone di lavaggio concentrato/ Tampón de lavado concentrado/ Tampão de Lavagem Concentrado/ wasbuffer, geconcentreerd/ Vaskebufferkoncentrat/ Vaskebufferkoncentrat/ tvättbuffertkoncentrat/ Bufor plukania koncentrat/ Mosópuffer koncentrátum/ Koncentrát vymývacieho pufru/ Концентрат на промивен буфер/ Pesurpuhvri kontsentrat/ Συμπύκνωμα ρυθμιστικού διαλύματος έκπλυσης/ Concentrat pentru tamponul de spălare/ Koncentrat izpiralnega pufru/ Pesuliuositiiviste |
| WASHBUF | Washing Buffer/ Waschpuffer/ Tampon de lavage/ Tampone di lavaggio/ Tampón de lavado/Tampão de Lavagem/ wasbuffer/ Vaskebuffer/ tvättbuffert/ Bufor plukania/ Mosópuffer/ Vymývací pufer/ Vymývací pufr/ Promivnen bufer/ Pesurpuhver/ Ρυθμιστικό διάλυμα έκπλυσης/ Tampon pentru spălare /Izpiralni pufer/ Pesuliuos |
| SUBST TMB | Substrate/ Substrat/ Substrat/ Substrato/ Substrato/ Substrato/ substraat/ Substrat/ Substrat/ Substrat/ Szubsztrátum/ Substrát/ Substrát/ Субстрат Substraat/ Υπόστρωμα/ Substrat/ Substrat/ Substraattiliuos |
| H₂SO₄ | Stopping Solution/ Stopplösung/ Stop Solution/ Soluzione di stop/ Stop Solución/ Solução Stop/ stopoplossing/ Stopopløsning/ Stopplösning/ Stop roztwór/ Megállító oldat/ Roztok na ukončenie/ Roztok pro ukončení/ Стопираш разтвор/ Stopp-lahus/ Διάλυμα διακοπής/ Soluție de oprire/ Stop raztopina/ Pysäytysliuos |
| TAPE | Cover Plate with sealing tape/ Platte abkleben/ Recouvrir la microplaque avec bande adhésive/ Coprire la piastra con nastro adesivo/ Cubrir la placa con una cinta adhesiva/ Cobrir a Placa com fita adesiva/ plaatje met tape afdekken/ Afdækningsplade med tape/ maskera platta/ Odkleić płytkę/ Tányér leragasztása/ Oblepit' podložku lepiacou páskou/ Olepiti podložku lepící páskou/ Плака с лента за запечатване/ Katta plaat isoleerikleepindiga/ Κολλήστε το πλακίδιο με κολλητική ταινία/ Aoperiți placa cu o bandă adezivă/ Prelepiti ploščo/ Peitã mikrotitrauslevy oheisella teipillä |
| MEASURE | Measure plate within 30 min at 450 nm (Referencefilter ≥590 nm)/ Ausmessung innerhalb von 30 min bei 450 nm (Referenzfilter ≥ 590 nm)/ Mesure lábsorbance en léspace de 30 min à 450 nm avec ≥590 nm longueur d'onde pour référence/ Misurazione entro 30 min. a 450 nm (filtro di riferimento ≥ 590 nm)/ Medición de la placa dentro de los siguientes 30 min a 450 nm (filtro de referencia ≥ 590nm)/ Medir a placa dentro de 30 min a 450 nm (Filtro de referência ≥ 590 nm)/ Binnen 30 minuten bij 450 nm meten (referentiefilter ≥ 590 nm)/ Mål plade i løbet af 30 min ved 450 nm (referencefilter ≥590 nm)/ Mät inom 30 min vid 450 nm (referensfilter ≥ 590 nm)/ Pomiar w ciągu 30 min przy 450 nm (filtr odniesienia ≥ 590 nm)/ Ki mérés 30 percen belül 450 nm-nél (referenciaszűrő ≥ 590 nm)/ Merať 30 minút pri 450 nm (Referenčných filtrov ≥590 nm)/ Měřit 30 minut při 450 nm (Referenční filtr ≥ 590 nm)/ Отчитане в рамките на 30 min при 450 nm (референтен филтър ≥ 590 nm)/ Mõõtmise 30 min jooksul 450 nm korral (võrdlusfilter ≥ 590 nm)/ Μέτρηση εντός 30 min στα 450 nm (φίλτρο αναφοράς ≥ 590 nm)/ Măsurare în decurs de 30 min la 450 nm (filtru de referință ≥ 590 nm)/ Izmerite ploščico v 30 min pri 450 nm (referenčni filter ≥590nm)/ Mittaa 30 minuutin aikana 450 nm:ssä (referenssi suodatin ≥ 590 nm) |
| Literatur | Literature/ Literatur/ Bibliographie/ Letterario/ Bibliografía/ Literatura documentação/ literatuur/ Litteratur/ litteratur/ Literatura/ Irodalom/ Literatúra/ Literatura/ Литература// Kirjandus/ Βιβλιογραφία/ Bibliografie/ literatura/ Lähdeluettelo |
| International Test description | International test description/ internationale Testanleitung/ description internationale de test/ Istruzioni per il test internazionali/ Descripción de ensayo internacional/ Descrição internacional do teste/ internationale testbeschrijving/ internationell testbeskrivning/ Opis testu międzynarodowego/ nemzetközi teszt-útmutató/ Medzinárodný návod k testu/ Mezinárodní návod k testu/ rahvusvaheline katse kirjeldus/ Διεθνείς οδηγίες για εργαστηριακές δοκιμές/ instructiuni internaționale pentru testare/ mednarodna navodila za preizkus/ Kansainvälinen käyttöohje |
| End | in all required wells/ in allen benötigten Vertiefungen/ dans tous les godets requis/ in tutti i pozzetti richiesti/ en todos los pozos requeridos/ em todos os tubos necessários/ in alle nodige putjes/ i alle nødvendige brønde/ i alla nödvändiga brunnar/ we wszystkich potrzebnych wgłębieniach/ minden szükséges forrásban/ vo všetkých potrebných miestach/ ve všech potřebných místech/ във всички необходими ямки/ kõigis vajalikes süvendites/ σε όλες τις απαραίτητες κοιλότητες/ în toate cavitățile necesare/ v vseh zahtevanih vdolbinah/ kaikkiin tarvittaviin mikrotitrauslevyn syvennyksiin |

TABLE OF CONTENTS

INSTRUCTIONS FOR USE

| | | |
|-----------|---|-----------|
| 1 | INTRODUCTION | 5 |
| 2 | INTENDED USE | 6 |
| 3 | ASSAY PRINCIPLE | 6 |
| 4 | WARNINGS AND PRECAUTIONS | 6 |
| 5 | SAMPLES | 7 |
| 6 | MATERIALS..... | 7 |
| 7 | TECHNICAL NOTES..... | 8 |
| 8 | ASSAY PROCEDURE..... | 9 |
| 9 | CALCULATION OF RESULTS | 10 |
| 10 | PERFORMANCE CHARACTERISTICS..... | 12 |
| 11 | LITERATURE / LITERATUR..... | 13 |
| 12 | INTERNATIONAL ASSAY DESCRIPTION..... | 14 |
| 13 | ASSAY PROCEDURE..... | 15 |

PACKAGE INSERT

| | |
|------------------------------|--|
| m/rIGFBP-3 ELISA E031 | 96 Bestimmungen |
| RUO | For Research Use Only |
| Principle of the test | Enzyme-linked Immunoassay |
| Duration (Incubation period) | 2.5 h |
| Antibodies | specific, goat anti-mouse/rat IGFBP-3 Antibody |
| Buffer | Ready for use and 20fold concentrate |
| Standard | 7 Single standards: (0.39 - 25 ng/mL), native Mouse IGFBP-3 |
| Assay Range | 0.09 – 12625 ng/mL |
| Control | 2 Control sera, lyophilized |
| Sample | Mouse- and Rat-Serum /Plasma Cell Culture Medium |
| Required sample volume | 10 µL |
| Sample dilution | 1:505 |
| Analytical Sensitivity | Ø 0.09 ng/mL |
| Intra- / Interassay Variance | Ø ≤10 % |

1 INTRODUCTION

Growth Hormone, Insulin-like Growth Factors and their binding proteins build up an endocrine system regulating not only longitudinal growth in humans but also influencing a broad variety of other physiological and pathophysiological processes like energy metabolism or tumor growth. Most effects of Growth Hormone (GH) are exerted by Insulin-like Growth Factors (IGF) mainly produced by the liver but also locally by specific tissues. The effects of IGF are also regulated, specific binding proteins (IGFBP 1-7) regulate bioavailability of IGF. After proteolytic cleavage of the binding proteins IGF is set free and able to bind to its receptor. The autophosphorylation of this tyrosine kinase receptor activates intra cellular signalling cascades. Some of these IGFBPs not only regulate the availability of IGF but also exert IGF-independent effects on cell physiology.

IGFBP-3 is the most abundant IGFBP in circulation and therefore of special relevance in regulation of IGF effects. This is reflected by the indicative value of serum IGFBP-3 concentration in diagnostics of growth disturbances. Regulation is effected e.g. through nourishing situation; Different diets for example affect the IGFBP-3 concentration (Bielohuby et al, 2010). IGFBP-3 has also been shown to be able to induce apoptosis, promote tumor growth and inhibit cellular migration and metastasis dependent on tissue and tumor stage.

Mouse / rat models for in vivo experiments are often used for studies of IGF-dependent and independent effects of IGFBP-3, particularly in the field of tumor research. For this purpose Mediagnost offers the E031 as a reliable and sensitive test system for the determination of IGFBP-3 in mouse and rat samples.

2 INTENDED USE

This enzyme immunoassay kit is suited for measuring IGFBP-3 in mouse and rat serum and plasma and in cell culture medium.

3 ASSAY PRINCIPLE

The Mediagnost m/rIGFBP-3 ELISA, E031 is a so-called sandwich-assay. It utilizes two different specific high affinity polyclonal antibodies for this protein. The IGFBP-3 in the samples binds quantitatively to the immobilized antibody. In the following step, the biotinylated antibody in turn binds IGFBP-3. After washing, a streptavidin-peroxidase-enzyme conjugate will be added, which will bind highly specific to the biotin of the antibody. Subsequently, the peroxidase catalyzes an enzymatic reaction resulting in a blue coloration. The intensity of the blue color depends on the IGFBP-3 content of the sample. The reaction is stopped by the addition of stop solution and color intensity is quantified by measuring the absorption.

4 WARNINGS AND PRECAUTIONS

For Research Use only. For Professional use only.

The Mediagnost kit is suitable only for in vitro and not for internal use in humans and animals. Follow strictly the test protocol. Use the valid version of the package insert provided with the kit. Be sure that everything has been understood. Mediagnost will not be held responsible for any loss or damage (except as required by statute) howsoever caused, arising out of noncompliance with the instructions provided.

Do not use obvious damaged or microbial contaminated or spilled material.

Caution: This kit contains material of human and/or animal origin. Therefore all components and patient's specimens should be treated as potentially infectious.

Appropriate precautions and good laboratory practices must be used in the storage, handling and disposal of the kit reagents. The disposal of the kit components must be made according to the local regulations.

Animal serum: mouse / rat in the following components: KS1, KS2

Reagents AK, EK, VP, WP

Contain as preservative **5-chloro-2-methyl-4-isothiazolin-3-one and 2-methyl-4-isothiazolin-3-one** (<0.015%)

| | |
|-----------|---|
| H317 | May cause an allergic skin reaction. |
| P280 | Wear protective gloves/ protective clothing/ eye protection/ face protection. |
| P272 | Contaminated work clothing should not be allowed out of the workplace. |
| P261 | Avoid breathing dust/ fume/ gas/ mist/ vapours/ spray. |
| P333+P313 | If skin irritation or rash occurs: Get medical advice/ attention. |
| P302+P352 | IF ON SKIN: Wash with plenty of soap and water. |
| P501 | Dispose of contents/ container in accordance with local/ regional/ national/ international regulations. |

Substrate Solution (S)

The TMB-Substrate (S) contains 3,3',5,5' Tetramethylbencidine (<0.05%)

| | |
|------------|---|
| H315 | Causes skin irritation. |
| H319 | Causes serious eye irritation. |
| H335 | May cause respiratory irritation. |
| P261 | Avoid breathing dust/ fume/ gas/ mist/ vapours/ spray. |
| P305+P351+ | IF IN EYES: Rinse cautiously with water for several minutes. |
| P338 | Remove contact lenses, if present and easy to do. Continue rinsing. |

Stopping Solution (SL)

The Stopping solution contains 0.2 M acid sulphur acid (H₂SO₄)

| | |
|------------|---|
| H290 | May be corrosive to metals. |
| H314 | Causes severe skin burns and eye damage. |
| P280 | Wear protective gloves/ protective clothing/ eye protection/ face protection. |
| P301+P330+ | IF SWALLOWED: rinse mouth. |
| P331 | Do NOT induce vomiting. |
| P305+P351+ | IF IN EYES: Rinse cautiously with water for several minutes. |
| P338 | Remove contact lenses, if present and easy to do. Continue rinsing. |
| P309+P310 | IF exposed or if you feel unwell: Immediately call a POISON CENTER or doctor/physician. |

4.1 General first aid procedures:

Skin contact: Wash affected area rinse immediately with plenty of water at least 15 minutes. Remove 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: After swallowing the product, if the affected person is conscious, rinse out the mouth with plenty of water: seek medical advice immediately.

5 SAMPLES

5.1 Sample type

Mouse and Rat Serum Plasma, In Heparin-Plasma samples the levels were found approx. 15% decreased. Further, cell culture medium was found to be suitable.

5.2 Specimen collection

Haemolytic reactions have to be avoided.

5.3 Requested sample volume: 10 µl serum.

5.4 Sample stability

- In firmly closable sample vials
- Storage at -20°C: min. 2 years
- Freeze/-thaw cycles: max. 3

It is recommended to store samples as soon as possible at least at 4°C. For any long time storage the sample has to be kept frozen at -20°C.

5.5 Sample dilution

Samples **must be diluted** prior to measurement. An extraction step is not required.

- Dilution: **1:505** with Dilution Buffer **VP**

We recommend a dilution in 2 steps:

Pipette **1 mL** Dilution Buffer **VP** in PE-/PP-Tubes (application of a multi-stepper is recommended in larger series), add **10 µL** Serum- or Plasma (dilution 1:101) and mix each tube immediately. Pipette **100 µL** of this dilution into another PE/PP vessel with **400 µL** of Dilution Buffer **VP** and mix immediately. This results in a final dilution of 1:505. After mixing, use 100 µL per assay in the assay within 1 hour of this solution


- After Mixing use 100 µL per assay in the assay within 1 hour of this solution.
- Where required, depending on the expected IGFBP-3-values, the dilution with **Dilution Buffer VP** can be higher or lower.
-

6 MATERIALS

6.1 Materials provided

The reagents listed below are sufficient for 96 wells including the standard curve.

| | | |
|------------|---|---------------------|
| MTP | Microtiter plate , ready for use, coated with goat anti-mouse IGFBP-3 Antibody, wells are separately breakable. | (8x12) wells |
| A-G | Standards , lyophilised (native Mouse-IGFBP-3), Concentrations are given on the vial labels and quality certificate. | 7 x 750 µL |
| KS1 | Control Serum 1 , lyophilised, (Mouse Serum), Concentration is given on the quality certificate . | 1x 250 µL |
| KS2 | Kontrollserum 2 , lyophilisiert, (Ratten Serum), Concentration is given on the quality certificate . | 1x 250 µL |
| AK | Antibody Conjugate , ready for use, Goat anti-mouse-IGFBP-3-Antibody, biotinylated . | 1 x 12 mL |
| EK | Enzyme Conjugate EK , contains HRP (Horseradish-Peroxidase)-labeled Streptavidin. | 1 x 12 mL |

| | | |
|---|--|-------------------|
| VP | Dilution Buffer , ready for use. Please shake before use. | 1 x 125 mL |
| WP | Washing Buffer WP , 20fold concentrated solution | 1 x 50 mL |
| S | Substrate S , ready for use, horseradish-peroxidase (HRP)- substrate, stabilised Tetramethylbencidine. | 1 x 12 mL |
| SL | Stopping Solution SL , ready for use, 0.2 M sulphuric acid. | 1 x 12 mL |
| - | Sealing Tape for covering the microtiter plate | 3 x |
|  | Instructions for use | 1 x |
| - | Quality Control Certificate (QC-Certificate) | 1 x |

6.2 Materials required, but not provided

- Distilled (Aqua destillata) or deionized water for dilution of the Washing Buffer **WP** (**A. dest.**), 950 mL.
- Graduated cylinder for diluting Washing Buffer (WP)
- Precision pipettes and multichannel pipettes with disposable plastic tips
- Polyethylene PE/Polypropylene PP tubes for dilution of samples
- Vortex-mixer
- Microtiter plate shaker (350 rpm)
- Microtiter plate washer (recommended)
- Micro plate reader ("ELISA-Reader") with filter for 450 and ≥ 590 nm
-

7 TECHNICAL NOTES

Storage Conditions

Store the kit at 2-8°C after receipt until its expiry date. The lyophilized reagents should be stored at -20 °C after reconstitution. Avoid repeated thawing and freezing.

Storage Life

The shelf life of the components **after initial opening** is warranted for **4 weeks**, store the unused strips and microtiter wells **airtight** together with the desiccant at 2-8°C in the clip-lock bag, use in the frame provided. The **reconstituted components** standards **A-G** and Control Sera **KS1 and KS2** must be stored at -20°C (max. 4 weeks). For further use, thaw quickly but gently (avoid temperature increase above room temperature and avoid excessive vortexing). Up to 3 of the freeze-thaw cycles did not influence the assay. The 1:20 diluted Washing Buffer **WP** is 4 weeks stable at 2-8°C

Preparation of reagents

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. Reagents with different lot numbers cannot be mixed.

Reconstitution

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

Dilution

After reconstitution dilute the Control Sera **KS1 and KS2** with the Dilution Buffer **VP** in the same ratio (1:505) as the sample. The required volume of Washing Buffer **WP** is prepared by 1:20 dilution of the provided 20-fold concentrate with Aqua dest.

Incubation

Incubation at room temperature means: Incubation at 20 - 25°C. The Substrate Solution **S**, stabilised H₂O₂-Tetramethylbencidine, is photosensitive—store and incubation in the dark.

Assay Procedure

When performing the assay, Blank, Standards **A-G**, Control Serum **KS** and the samples should be pipette as fast as possible (e.g. <15 minutes). To avoid distortions due to differences in incubation times, Antibody Conjugate **AK**, Enzyme Conjugate **EK** as well as the succeeding Substrate Solution **S** should be added to the plate in the same order and in the same time interval as the samples. Stopping Solution **SL** should be added to the plate in the same order as Substrate Solution **S**.

All determinations (Blank, Standards **A-G**, Control Serum **KS** and samples) should be assayed in duplicate. For optimal results, accurate pipetting and adherence to the protocol are recommended.

Shaking

The incubation steps should be performed at mean rotation frequency of a particularly suitable microtiter plate shaker. We are recommending 350 rpm. Due to certain technical differences deviations may occur, in case the rotation frequency must be 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.

Washing

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 **WP** 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 microtiter** 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 dynamical swinging out the microtiter 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.

8 ASSAY PROCEDURE

| Reagent preparation | | Reconstitution | Dilution |
|--|--|--|---|
| A-G | Standards | in 750 µL Dilution Buffer VP | - |
| KS1 | Control Serum 1 | in 250 µL Dilution Buffer VP | 1:505 with Dilution Buffer VP |
| KS2 | Control Serum 2 | in 250 µL Dilution Buffer VP | 1:505 with Dilution Buffer VP |
| WP | Waschpuffer | - | 1:20 with Aqua dest. |
| Dilute Samples with Dilution Buffer VP 1:505 | | | |
| Before assay procedure bring all reagents to room temperature (20°C- 25°C) | | | |
| Assay Procedure in Double Determination: | | | |
| Pipette | Reagents | | Position |
| 100 µL | Dilution Buffer VP (Blank) | | A1/A2 |
| 100 µL | Standard A (0.39 ng/mL) | | B1/B2 |
| 100 µL | Standard B (0.78 ng/mL) | | C1/C2 |
| 100 µL | Standard C (1.56 ng/mL) | | D1/D2 |
| 100 µL | Standard D (3.13 ng/mL) | | E1/E2 |
| 100 µL | Standard E (6.25 ng/mL) | | F1/F2 |
| 100 µL | Standard F (12.5 ng/mL) | | G1/G2 |
| 100 µL | Standard G (25 ng/mL) | | H1/H2 |
| 100 µL | Control Serum KS 1 | (1:505 diluted) | A3/A4 |
| 100 µL | Control Serum KS 2 | (1:505 diluted) | B3/B4 |
| 100 µL | Sample | (1:505 diluted) | In the rest of the wells according to the requirements. |
| Cover the wells with the sealing tape. | | | |
| Incubation: 1 h at 20-25°C, 350 rpm | | | |
| 5x 300 µL | Aspirate the contents of the wells and wash 5x with 300 µL each WP/well | | Each well |
| 100 µL | Antibody Conjugat AK | | Each well |
| Cover the wells with the sealing tape | | | |

| Incubation: 1 h at 20-25°C, 350 rpm | | |
|--|--|-----------|
| 5x 300 µL | Aspirate the contents of the wells and wash 5x with 300 µL each WP/well | Each well |
| 100 µL | Enzyme Conjugate EK | Each well |
| Cover the wells with the sealing tape | | |
| Incubation: 15 min at 20-25°C, 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 |
| Substrat S Incubation: 15 Minutes in the dark at RT | | |
| 100 µL | Stop Solution SL | Each well |
| Measure the absorbance within 30 min at 450 nm (≥ 590 nm Reference) | | |

9 CALCULATION OF RESULTS

For the evaluation of the assay it is required that the absorbance values of the blank should be below 0.25, and the absorbance of standard G should be above 1.00.

Samples, which yield higher absorbance values than **Standard G**, are beyond the standard curve, for reliable determinations such samples should be retested at a higher dilution.

9.1 Establishing the standard curve

The standards provided contain the following concentration of **mIGFBP-3**:

| Standard | A | B | C | D | E | F | G |
|----------|------|------|------|------|------|------|----|
| ng/mL | 0.39 | 0.78 | 1.56 | 3.13 | 6.25 | 12.5 | 25 |

- 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 **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 5.760 ng/mL, Sample was 1:505 diluted: $5.760 \times 505 = 2909$ ng/mL, or 2.909 µg/mL equal to 2.909 mg/L)

9.2 Example of a typical standard curve

| Standard | Blank | A | B | C | D | E | F | G |
|-----------------|-------|-------|-------|-------|-------|-------|-------|-------|
| ng/mL | 0 | 0.39 | 0.78 | 1.56 | 3.13 | 6.25 | 12.5 | 25 |
| OD (450-620 nm) | 0.00 | 0.048 | 0.101 | 0.202 | 0.412 | 0.815 | 1.484 | 2.438 |

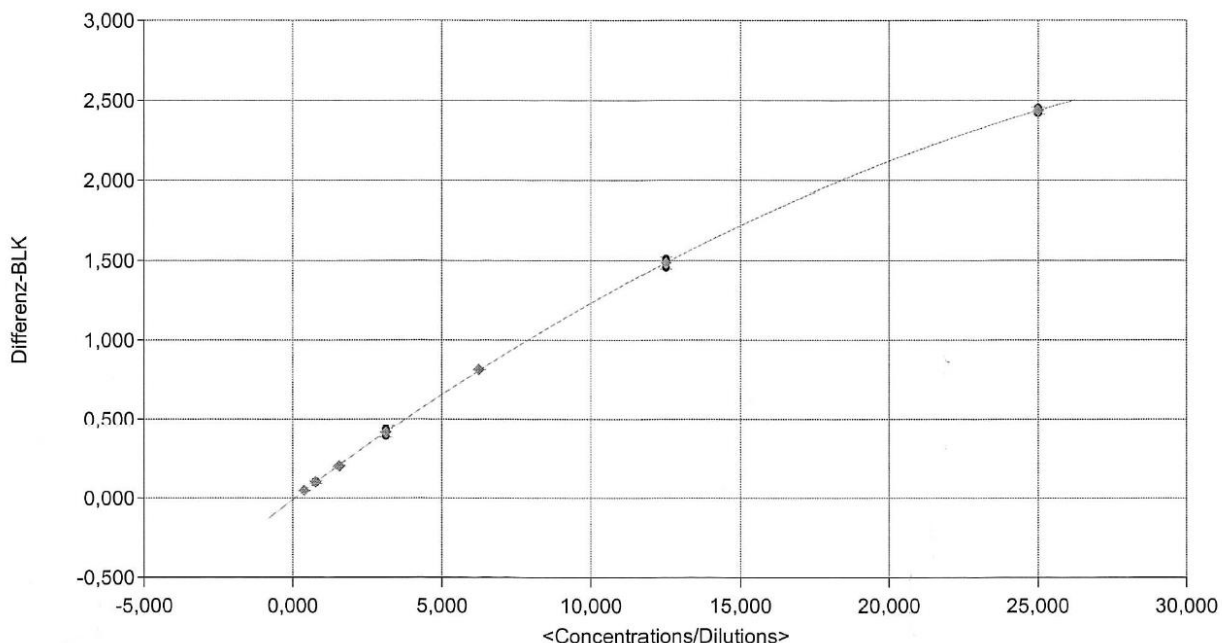


Fig. 1: Exemplary Standard Curve with a polynomial 3rd degree as curve fit.

The exemplary shown standard curve in Fig.1 **cannot be used** for calculation of your test results. You have to establish a standard curve for each test you conduct!

Exemplary calculation of the IGFBP-3 concentration of a diluted sample:

OD 450 nm

Measured extinction (mean value) of your sample 0.749

Measured extinction of the blank (mean value) 0.000

Your **measurement program** will calculate the IGFBP-3 concentration of the sample automatically by using the difference of sample and blank for the calculation. You only have to determine the most suitable curve fit (here: polynomial 3rd degree).

In this exemplary case the following equation is solved by the program to calculate the IGFBP-3 concentration in the sample:

$$0.749 = 3,62^{-6} \times X^3 + -0.00188 \times X^2 + 0.143 \times X - 0.011$$
$$5.760 = X$$

Multiplication by dilution factor (1:505) gives the IGFBP-3 concentration of the sample with 2909 ng/mL

10 PERFORMANCE CHARACTERISTICS

10.1 Calibration

The Mediagnost E031 was recalibrated on a highly purified eukaryotic expressed recombinant rat IGFBP-3 preparation. The previous calibration on a recombinant mouse IGFBP-3 preparation of lower purity and with a partly different amino acid sequence resulted approx. by a factor of 10 lower values. If desired for comparison purposes, a conversion from or to old values can be carried out by a factor of 10. Measured values with **previous calibration**, e.g. 300 ng/mL correspond to a **new calibration** of 3000 ng/mL. It is also possible to calculate the previous calibration by dividing by 10.

10.2 Analytical Sensitivity

The **analytical Sensitivity** was assessed by 21-fold determination of the blank and calculating the theoretical concentration of the blank +2SD. The analytical sensitivity of the E031 is **0.09 ng/mL**.

10.3 Precision

The **Inter-** and **Intra-Assay** variation coefficients were on average $\leq 10\%$. Exemplary determinations are shown in table 1 and table 2.

Tabelle 1 Inter-Assay-Variation (n=26 or 15)

| | Mean Value (ng/mL) | Standard Deviation (ng/mL) | VC(%) |
|----------|--------------------|----------------------------|-------|
| Sample 1 | 4836 | 384 | 7,94 |
| Sample 2 | 2625 | 269 | 9,43 |

Tabelle 2 Intra-Assay-Variation (n=13)

| | Mean Value (ng/mL) | Standard Deviation (ng/mL) | VC (%) |
|----------|--------------------|----------------------------|--------|
| Sample 1 | 3286 | 124 | 3,76 |
| Sample 2 | 1529 | 123 | 8,02 |

10.4 Linearity

Table 3: Linearity (results of 2 different mouse sera)

| Dilution: | Sample 1 (recalculated, ng/mL) | Sample 2 (recalculated, ng/mL) |
|--------------------|-----------------------------------|-----------------------------------|
| 1:100 | 3518 | 3676 |
| 1:200 | 3691 | 4145 |
| 1:400 | 3845 | 4234 |
| 1:800 | 3813 | 4110 |
| 1:1600 | 3792 | 4219 |
| 1:3200 | 3861 | 4557 |
| AV/ SD/ VC% | 3753/ 129/ 3.46 | 4157/ 284/ 6.83 |

AV = Average Value, SD = Standard Deviation; VC = Coefficient of Variation

10.5 Species Cross-Reactivity

Serum of the cited species were used as diluted samples in this assay system.

No cross reactivity was detected for:

Rabbit, Cat, Chicken, Guinea pig, Goat, Sheep, Horse, Donkey, Pig, Dog, Bovine.

Cross reactivity with recombinant human eukaryotic expressed IGFBP-3 (1 $\mu\text{g/mL}$): 0.06%

10.6 Recovery in Cell Culture Medium

The recovery of recombinant mouse IGFBP-3 in **cell culture medium** DMEM was found to be 89.4%, and, in DMEM incl. 5% FCS 92.6%. Therefore, cell culture medium seems to be suitable as sample matrix.

11 LITERATURE / LITERATUR

Silha JV, Murphy LJ: Insulin-like growth factor binding proteins in development. *Adv Exp Med Biol.* 2005;567:55-89

Yakar S, Sun H, Zhao H, Pennisi P, Toyoshima Y, Setser J, Stannard B, Scavo L, Leroith D: Metabolic effects of IGF-I deficiency: lessons from mouse models. *Pediatr Endocrinol Rev.* 2005 Sep;3(1):11-19

Silha JV, Murphy LJ: Insights from insulin-like growth factor binding protein transgenic mice. *Endocrinology.* 2002 Oct;143(10):3711-4.

Firth SM, Baxter RC: Cellular actions of the insulin-like growth factor binding proteins. *Endocr Rev.* 2002 Dec;23(6):824-854

Schneider MR, Lahm H, Wu M, Hoefflich A, Wolf E: Transgenic mouse models for studying the functions of insulin-like growth factor-binding proteins. *FASEB J.* 2000 Apr;14(5):629-40

Bielohuby M, Matsuura M, Herbach N, Kienzle E, Slawik M, Hoefflich A, Bidlingmaier M. Short Term Exposure to Low-Carbohydrate / High Fat Diets Induces Low Bone Mineral Density and Reduces Bone Formation in Rats. *Journal of Bone and Mineral Research*, Vol. 25, No. 2, February 2010, pp 275–284

INTERNATIONAL ASSAY DESCRIPTION

| | | | | |
|--|-----------------------------------|-------------------|--------|--------------------|
| A-G | STD | Rec in 750 µL | BUF VP | 100 µL |
| KS1 | Control | Rec in 250 µL | BUF VP | 100 µL |
| KS2 | Control | Rec in 250 µL | BUF VP | 100 µL |
| WP | WASHBUF 20x | - | | 1:20 DILU A. dest. |
| SPE + Control 1:505 DILU BUF VP ↔ ⌚ max. 1 h | | | | 100 µL |
| °C 20-25°C | | | | |
| 100 µL | BUF VP | | | A1/2 |
| 100 µL | STD A | (0.39 ng/mL) | | B1/2 |
| 100 µL | STD B | (0.78 ng/mL) | | C1/2 |
| 100 µL | STD C | (1.56 ng/mL) | | D1/2 |
| 100 µL | STD D | (3.13 ng/mL) | | E1/2 |
| 100 µL | STD E | (6.25 ng/mL) | | F1/2 |
| 100 µL | STD F | (12.5 ng/mL) | | G1/2 |
| 100 µL | STD G | (25 ng/mL) | | H1/2 |
| 100 µL | CONTROL KS1 | 1:505 DILU BUF VP | | A3/4 |
| 100 µL | CONTROL KS2 | 1:505 DILU BUF VP | | B3/4 |
| 100 µL | SPE 1:505 | DILU BUF VP | | |
| TAPE | | | | |
| ⌚ 1 h °C 20-25 ↔ 350 rpm | | | | |
| 5x 300 µL | 5x WASHBUF WP | | | |
| 100 µL | Ab AK | | | |
| TAPE | | | | |
| ⌚ 1 h °C 20-25 ↔ 350 rpm | | | | |
| 5x 300 µL | 5x WASHBUF WP | | | |
| 100 µL | CONJ EK | | | |
| TAPE | | | | |
| ⌚ 15 min °C 20-25 ↔ 350 rpm | | | | |
| 5x 300 µL | 5x WASHBUF WP | | | |
| 100 µL | SUBST TMB S | | | |
| ⌚ 15 min °C 20-25 ☀ | | | | |
| 100 µL | H ₂ SO ₄ SL | | | |
| MEASURE | | | | |

12 ASSAY PROCEDURE

| Reagent preparation | | Reconstitution | Dilution |
|--|--|--|---|
| A-G | Standards | in 750 µL Dilution Buffer VP | - |
| KS1 | Control Serum 1 | in 250 µL Dilution Buffer VP | 1:505 with Dilution Buffer VP |
| KS2 | Control Serum 2 | in 250 µL Dilution Buffer VP | 1:505 with Dilution Buffer VP |
| WP | Waschpuffer | - | 1:20 with Aqua dest. |
| Dilute Samples with Dilution Buffer VP 1:505 | | | |
| Before assay procedure bring all reagents to room temperature (20°C- 25°C) | | | |
| Assay Procedure in Double Determination: | | | |
| Pipette | Reagents | | Position |
| 100 µL | Dilution Buffer VP (Blank) | | A1/A2 |
| 100 µL | Standard A (0.39 ng/mL) | | B1/B2 |
| 100 µL | Standard B (0.78 ng/mL) | | C1/C2 |
| 100 µL | Standard C (1.56 ng/mL) | | D1/D2 |
| 100 µL | Standard D (3.13 ng/mL) | | E1/E2 |
| 100 µL | Standard E (6.25 ng/mL) | | F1/F2 |
| 100 µL | Standard F (12.5 ng/mL) | | G1/G2 |
| 100 µL | Standard G (25 ng/mL) | | H1/H2 |
| 100 µL | Control Serum KS 1 | (1:505 diluted) | A3/A4 |
| 100 µL | Control Serum KS 2 | (1:505 diluted) | B3/B4 |
| 100 µL | Sample | (1:505 diluted) | In the rest of the wells according to the requirements. |
| Cover the wells with the sealing tape. | | | |
| Incubation: 1 h at 20-25°C, 350 rpm | | | |
| 5x 300 µL | Aspirate the contents of the wells and wash 5x with 300 µL each WP/well | | Each well |
| 100 µL | Antibody Conjugat AK | | Each well |
| Cover the wells with the sealing tape | | | |
| Incubation: 1 h at 20-25°C, 350 rpm | | | |
| 5x 300 µL | Aspirate the contents of the wells and wash 5x with 300 µL each WP/well | | Each well |
| 100 µL | Enzyme Conjugate EK | | Each well |
| Cover the wells with the sealing tape | | | |
| Incubation: 15 min at 20-25°C, 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 |
| Substrat S Incubation: 15 Minutes in the dark at RT | | | |
| 100 µL | Stop Solution SL | | Each well |
| Measure the absorbance within 30 min at 450 nm (≥590 nm Reference) | | | |