

**Manual**

# **IDK® DAO ELISA**

***For the in vitro determination of DAO  
in serum and dried blood spots***

Valid from 2025-04-22

**REF K 8500**

**REF K 8500.20**

$\Sigma$  96

$\Sigma$  20 x 96



**Immundiagnostik AG**, Stubenwald-Allee 8a, 64625 Bensheim, Germany

Tel.: +49 6251 70190-0

Fax: + 49 6251 70190-363

e.mail: [info@immundiagnostik.com](mailto:info@immundiagnostik.com)

[www.immundiagnostik.com](http://www.immundiagnostik.com)

# Table of Contents

|   |           |
|---|-----------|
| <b>1. INTENDED USE</b>                                  | <b>21</b> |
| <b>2. INTRODUCTION</b>                                  | <b>21</b> |
| <b>3. MATERIAL SUPPLIED</b>                             | <b>22</b> |
| <b>4. MATERIAL REQUIRED BUT NOT SUPPLIED</b>            | <b>23</b> |
| <b>5. STORAGE AND PREPARATION OF REAGENTS</b>           | <b>23</b> |
| <b>6. STORAGE AND PREPARATION OF SAMPLES</b>            | <b>24</b> |
| <i>Preanalytic handling</i>                             | 24        |
| <i>Serum</i>  | 24        |
| <i>Dried blood spots</i>                                | 24        |
| <b>7. ASSAY PROCEDURE</b>                               | <b>25</b> |
| <i>Principle of the test</i>                            | 25        |
| <i>Test procedure</i>                                   | 25        |
| <b>8. RESULTS</b>                                       | <b>27</b> |
| <b>9. LIMITATIONS</b>                                   | <b>28</b> |
| <b>10. QUALITY CONTROL</b>                              | <b>28</b> |
| <i>Reference range</i>                                  | 28        |
| <i>Heparin treatment</i>                                | 29        |
| <b>11. PERFORMANCE CHARACTERISTICS</b>                  | <b>30</b> |
| <i>Accuracy – Precision</i>                             | 30        |
| <i>Accuracy – Trueness</i>                              | 31        |
| <i>Linearity</i>  | 32        |
| <i>Analytical sensitivity</i>                           | 32        |
| <b>12. PRECAUTIONS</b>                                  | <b>33</b> |
| <b>13. TECHNICAL HINTS</b>                              | <b>33</b> |
| <b>14. GENERAL NOTES ON THE TEST AND TEST PROCEDURE</b> | <b>34</b> |
| <b>15. REFERENCES</b>                                   | <b>34</b> |

## 1. INTENDED USE

This Immundiagnostik AG assay is an enzyme immunoassay intended for the quantitative determination of diamine oxidase (DAO) in serum and dried blood spots.

For *in vitro* diagnostic use only.

## 2. INTRODUCTION

Diamine oxidase (DAO) is a body's own enzyme that metabolises histamine. Although DAO is found practically in the whole body, the most important site of its action is the intestine. The enzymatic activity of DAO determines the histamine degradation speed. In the case of DAO deficiency or inhibition, incorporated or endogenous histamine cannot be degraded quickly enough, and the symptoms of histamine intolerance are presented. Millions of people suffer from gastrointestinal problems, migraine, irritations of nasal mucosa and other allergy-like symptoms after consumption of certain nutrients. Too much histamine in the body can be the reason for this wide range of symptoms.

The determination of DAO serum concentration (K 8500) combined with the determination of DAO activity (K 8220 DAO REA) is a suitable marker for the differential diagnosis of histamine intolerance and associated symptoms.

Our IDK® DAO ELISA kit is intended for determination of the diamine oxidase (DAO) concentration in serum.

### Indications

- Frequent headaches or migraine
- Snuffles after consumption of histamine-containing nutrients
- Tissue oedema
- Eyelid turgor
- Skin redness
- Limb aches
- Gastrointestinal discomfort
- Monitoring of a histamine free diet

### 3. MATERIAL SUPPLIED

| Cat. No      | Label      | Kit Components  | Quantity     |                   |
|--------------|------------|---|--------------|-------------------|
|              |            |   | K 8500       | K 8500.20         |
| K 8500       | PLATE      | Microtiter plate, pre-coated                              | 12 x 8 wells | 20 x 12 x 8 wells |
| K 0001.C.100 | WASH-BUF   | Wash buffer concentrate, 10x                              | 2 x 100 ml   | 40 x 100 ml       |
| K 8500       | STD        | Standards, lyophilised (10; 3.3; 1.1; 0.37; 0 U/ml)*      | 4 x 5 vials  | 25 x 5 vials      |
| K 8500       | CTRL1      | Control, lyophilised (see specification for range)        | 4 x 1 vial   | 25 x 1 vial       |
| K 8500       | CTRL2      | Control, lyophilised (see specification for range)        | 4 x 1 vial   | 25 x 1 vial       |
| K 8500       | AB         | Detection antibody concentrate, biotinylated              | 1 x 200 µl   | 20 x 200 µl       |
| K 8500       | CONJ       | Conjugate concentrate, peroxidase-labelled (streptavidin) | 1 x 200 µl   | 20 x 200 µl       |
| K 8500       | ABBUF      | Dilution buffer for AB and CONJ, ready-to-use             | 1 x 50 ml    | 20 x 50 ml        |
| K 8500       | SAMPLE-BUF | Sample dilution buffer, ready-to-use                      | 1 x 50 ml    | 20 x 50 ml        |
| K 0002.15    | SUB        | Substrate (tetramethylbenzidine), ready-to-use            | 1 x 15 ml    | 20 x 15 ml        |
| K 0003.15    | STOP       | Stop solution, ready-to-use                               | 1 x 15 ml    | 20 x 15 ml        |

For reorders of single components, use the catalogue number followed by the label as product number.

**\*IMPORTANT: For information on the measuring range of the product, see chapters 8 and 9.**

#### 4. MATERIAL REQUIRED BUT NOT SUPPLIED

- Ultrapure water\*
- Calibrated precision pipettors and 10–1000 µl single-use tips
- Standard laboratory reaction vessels 1.5 ml (single-use)
- Standard laboratory reaction vessel (15 ml) (single-use)
- Foil to cover the microtiter plate
- Centrifuge, 3000 g
- Multi-channel pipets or repeater pipets
- Vortex
- Microtiter plate thermoshaker at 37 °C (for example model Shake ID2 available at Immundiagnostik AG)
- Microtiter plate reader (required filters see chapter 7)

\* Immundiagnostik AG recommends the use of ultrapure water (water type 1; ISO 3696), which is free of undissolved and colloidal ions and organic molecules (free of particles > 0.2 µm) with an electrical conductivity of 0.055 µS/cm at 25 °C (≥ 18.2 MΩ cm).

#### 5. STORAGE AND PREPARATION OF REAGENTS

- To run the assay more than once, ensure that reagents are stored at the conditions stated on the label. **Prepare only the appropriate amount necessary for each run.** The kit can be used up to 4 times within the expiry date stated on the label.
- Reagents with a volume less than **100 µl** should be centrifuged before use to avoid loss of volume.
- **Preparation of the wash buffer:** The **wash buffer concentrate (WASHBUF)** has to be diluted with ultrapure water **1:10** before use (100 ml WASHBUF + 900 ml ultrapure water), mix well. Crystals could occur due to high salt concentration in the concentrate. Before dilution, the crystals have to be redissolved at room temperature or in a water bath at 37 °C. The **WASHBUF** can be used until the expiry date stated on the label when stored at **2–8 °C**. **Wash buffer** (1:10 diluted WASHBUF) can be stored in a closed flask at **2–8 °C for 1 month**.
- The **lyophilised standards (STD)** and **controls (CTRL)** are stable at **2–8 °C** until the expiry date stated on the label. Before use, the STD and CTRL have to be reconstituted with **500 µl sample dilution buffer (SAMPLEBUF)** and mixed by gentle inversion to ensure complete reconstitution. Allow the vial content to dissolve for 10 minutes and then mix thoroughly. **Standards and controls** (reconstituted STD and CTRL) **are not stable and cannot be stored**.

- **Preparation of the conjugate and the detection antibody:** Before use, the **conjugate concentrate (CONJ)** and the **detection antibody concentrate (AB)** have to be diluted **1:101** in **dilution buffer** (100 µl CONJ + 10 ml ABBUF), (100 µl AB + 10 ml ABBUF). The CONJ and the AB are stable at **2–8 °C** until the expiry date stated on the label. **Conjugate** (1:101 diluted CONJ) and **detection antibody** (1:101 diluted AB) **are not stable and cannot be stored.**
- All other test reagents are ready-to-use. Test reagents are stable until the expiry date (see label) when stored at **2–8 °C**.

## 6. STORAGE AND PREPARATION OF SAMPLES

### *Preanalytic handling*

**Lipemic or hemolytic samples may give erroneous results and should not be used for analysis.**

### *Serum*

#### **Sample storage**

The samples can be stored for **6 months at -20 °C**. Avoid repeated freezing and thawing. The samples are stable at **room temperature** for up to **4 days** and at **2–8 °C** for up to **9 days**.

#### **Sample preparation**

Serum samples must be diluted **1:5** before performing the assay,

e.g. **50 µl** sample + **200 µl** sample dilution buffer (SAMPLEBUF), mix well.

**100 µl** of the dilution are used in the test per well.

### *Dried blood spots*

#### **Collection and storage of dried blood spots**

**50 µl whole blood** dripped on a dried sample carrier cleared by Immundiagnostik AG are suitable as sample material after complete drying. We recommend DrySpot-ID (catalogue no DZ9020ID or DZ9021ID) as dried blood spot carrier.

## Preparation of the dried blood samples

|    |  |
|----|--|
| 1. | Label 1,5- ml polypropylene tubes..  |
| 2. | Remove filter from sampling device (always wear gloves since the sample is potentially infectious).                                      |
| 3. | Put filter in a labelled tube.   |
| 4. | Add <b>400 µl</b> sample dilution buffer (SAMPLEBUF) per sample, allow sample to stand for <b>20 min</b> at room temperature (15–30 °C). |
| 5. | Vortex for <b>10 sec</b> . The filter will decolourise.  |
| 6. | Centrifuge the samples for <b>5 min</b> at <b>3000 g</b> to remove residual filter pieces.   |

## 7. ASSAY PROCEDURE

### *Principle of the test*

This ELISA is designed for the quantitative determination of DAO in serum and dried blood spots. The assay utilises the sandwich technique with two polyclonal antibodies against recombinant DAO.

Standards, controls and prepared samples which are assayed for DAO are added into the wells of a micro plate coated with polyclonal rabbit anti-DAO antibody. During the first incubation step, DAO is bound by the immobilised primary antibody. Then a biotinylated polyclonal anti-DAO antibody is added into each microtiter well. In the next step, the streptavidin peroxidase conjugate is added and a “sandwich” of 1st antibody – DAO – biotinylated antibody – streptavidin peroxidase conjugate is formed. Tetramethylbenzidine (TMB) is used as peroxidase substrate. Finally, an acidic stop solution is added to terminate the reaction. The colour changes from blue to yellow. The intensity of the yellow colour is directly proportional to the concentration of DAO. A dose response curve of the absorbance unit (optical density, OD at 450 nm) vs. concentration is generated, using the values obtained from the standard. DAO, present in the patient samples, is determined directly from this curve.

### *Test procedure*

Bring all **reagents and samples to room temperature** (15–30 °C) and mix well.

Mark the positions of standards/controls/samples on a protocol sheet.

Take as many microtiter strips as needed from kit. Store unused strips together with the desiccant bag in the closed aluminium packaging at 2–8 °C. Strips are stable until expiry date stated on the label.

For automated ELISA processors, the given protocol may need to be adjusted according to the specific features of the respective automated platform. For further details please contact your supplier or Immundiagnostik AG.

We recommend to carry out the tests in duplicate.

|     |   |
|-----|---|
| 1.  | <b>Before use</b> , wash the wells <b>5 times</b> with <b>250 µl wash buffer</b> . After the final washing step, remove residual wash buffer by firmly tapping the plate on absorbent paper.        |
| 2.  | Add each <b>100 µl standards/controls/prepared samples</b> into the respective wells.   |
| 3.  | Cover the strips and incubate for <b>2 hours</b> at <b>37 °C</b> on a <b>horizontal shaker*</b> .   |
| 4.  | Discard the content of each well and wash <b>5 times</b> with <b>250 µl wash buffer</b> . After the final washing step, remove residual wash buffer by firmly tapping the plate on absorbent paper. |
| 5.  | Add <b>100 µl detection antibody</b> (diluted AB) into each well, mix gently.   |
| 6.  | Cover the strips and incubate for <b>1 hour</b> at <b>37 °C</b> on a <b>horizontal shaker*</b> .  |
| 7.  | Discard the contents of each well and wash <b>5 times</b> with <b>250 µl wash buffer</b> . After the final washing step, the inverted microtiter plate should be firmly tapped on absorbent paper.  |
| 8.  | Add <b>100 µl conjugate</b> (diluted CONJ) into each well.  |
| 9.  | Cover the strips and incubate for <b>1 hour</b> at <b>37 °C</b> on a <b>horizontal shaker*..</b>  |
| 10. | Discard the contents of each well and wash <b>5 times</b> with <b>250 µl wash buffer</b> . After the final washing step, the inverted microtiter plate should be firmly tapped on absorbent paper.  |
| 11. | Add <b>100 µl substrate</b> (SUB) into each well.   |
| 12. | Incubate for <b>10–20 min**</b> at room temperature (15–30 °C) in the <b>dark</b> .   |
| 13. | Add <b>100 µl stop solution</b> (STOP) into each well and mix well.   |



|     |   |
|-----|---|
| 14. | Determine <b>absorption immediately</b> with an ELISA reader at <b>450 nm</b> against 620 nm (or 690 nm) as a reference. If no reference wavelength is available, read only at 450 nm. If the extinction of the highest standard exceeds the range of the photometer, absorption must be measured immediately at <b>405 nm</b> against 620 nm as a reference. |
|-----|---|

- \* We recommend shaking the strips at 700 rpm with an orbit of 2 mm.
- \*\* The intensity of the colour change is temperature sensitive. We recommend observing the colour change and stopping the reaction upon good differentiation.

8. RESULTS

The following algorithms can be used alternatively to calculate the results. We recommend using the “4 parameter algorithm”.

1. 4 parameter algorithm

It is recommended to use a linear ordinate for the optical density and a logarithmic abscissa for the concentration. When using a logarithmic abscissa, the zero standard must be specified with a value less than 1 (e.g. 0.001).

2. Point-to-point calculation

We recommend a linear ordinate for the optical density and a linear abscissa for the concentration.

3. Spline algorithm

We recommend a linear ordinate for the optical density and a linear abscissa for the concentration.

The plausibility of the duplicate values should be examined before the automatic evaluation of the results. If this option is not available with the programme used, the duplicate values should be evaluated manually.

Serum

The obtained results have to be multiplied by the **dilution factor of 5** to get the actual concentrations.

Dried blood spots

The obtained results have to be multiplied by the **factor of 6** to get the actual concentrations.

In case **another dilution factor** has been used, multiply the obtained result by the dilution factor used.

## 9. LIMITATIONS

Samples with concentrations above the measurement range (see definition below) can be further diluted and re-assayed. Please consider this higher dilution when calculating the results.

Samples with concentrations lower than the measurement range (see definition below) cannot be clearly quantified.

The upper limit of the measurement range can be calculated as:

*highest concentration of the standard curve × sample dilution factor to be used*

The lower limit of the measurement range can be calculated as:

*LoB × sample dilution factor to be used*

LoB see chapter "Performance Characteristics".

## 10. QUALITY CONTROL

Immundiagnostik AG recommends the use of external controls for internal quality control, if possible.

Control samples should be analysed with each run. Results, generated from the analysis of control samples, should be evaluated for acceptability using appropriate statistical methods. The results for the patient samples may not be valid if within the same assay one or more values of the quality control sample are outside the acceptable limits.

### *Reference range*

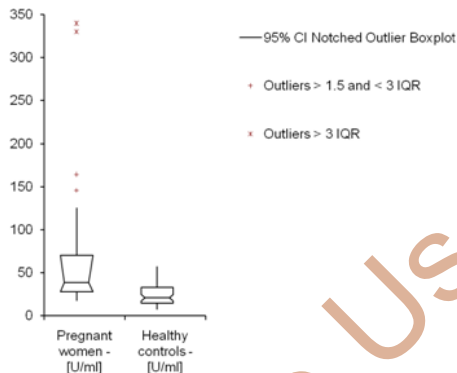
- |              |  |
|--------------|--|
| < 3 U/ml:    | high incidence for HIT (Histamine intolerance) |
| 3 - 10 U/ml: | HIT probable                                   |
| > 10 U/ml:   | low incidence for HIT                          |

Conversion factor: 1 U/ml = 1.25 ng/ml

We recommend each laboratory to establish its own reference range.

**Comparison “Pregnant women” and “Healthy controls”**

For the clinical evaluation of this assay we have analysed samples from pregnant women and apparently healthy controls. The IDK® DAO ELISA detects, as required and expected, higher values in pregnant women than in healthy controls.



**Heparin treatment**

Furthermore, DAO levels in healthy study participants increased sharply within 30 minutes of heparin administration. It has been documented in scientific literature that DAO levels rise after heparin administration.

**Treatment outcome before and after heparin administration**

IDK® DAO ELISA [U/ml]

|           | BEFORE administration | 30 min AFTER | 60 min AFTER |
|-----------|-----------------------|--------------|--------------|
| Patient 1 | 76                    | 219          | -            |
| Patient 2 | 55                    | 152          | -            |
| Patient 3 | 2.5                   | 277          | 622          |
| Patient 4 | 18.9                  | 621          | 555          |

Since the IDK® DAO ELISA determines DAO concentration while the conventional histamine intolerance tests with putrescine or histamine as substrate, e.g. DAO REA, K 8220.2, determine DAO activity, the correlation coefficient must not necessarily be  $r > 0.8$ . This can be explained by the fact that the activity does not depend on the number of molecules alone, but also on cofactors such as vitamin C, vitamin B6, copper or manganese ions in vitro and in vivo. For the diagnosis of histamine intolerance via DAO activity test we therefore recommend to determine the above mentioned cofactors as well. The problem may not be a low DAO level, but a cofactor deficiency.

The symptoms of histamine intolerance can be caused by low DAO activity because the above-mentioned cofactors are not sufficiently available. By quantitating the cofactors it can be determined which one needs to be supplemented.

### Medication effects

In addition, histamine intolerance symptoms may be due to low DAO activity caused by medication such as:

|                                |   |
|--------------------------------|---|
| Muscle relaxants               | Pancuronium, alcuronium, D-tubocurarine   |
| Narcotics                      | Thiopental  |
| Analgetics                     | Morphine, pethidine, nonsteroidal anti-inflammatory drugs, acetylsalicylic acid, metamizole |
| Local anesthetics              | Prilocaine  |
| Antihypotonics                 | Dobutamine  |
| Antihypertensive drugs         | Verapamil, alprenolol, dihydralazine  |
| Antiarrhythmics                | Propafenone   |
| Diuretics                      | Amiloride   |
| Drugs influencing gut motility | Metoclopramide  |
| Antibiotics                    | Cefuroxime, cefotiam, isoniazid, pentamidin, clavulanic acid, chloroquine                   |
| Mucolytics                     | Acetylcysteine, ambroxol  |
| Broncholytics                  | Aminophylline   |
| H2-receptor antagonists        | Cimetidine  |
| Cytostatics                    | Cyclophosphamide  |
| Antidepressants                | Amitriptyline   |

If you are taking such medication, you may want to discuss with your physician alternative medication in order to relieve your symptoms.

## 11. PERFORMANCE CHARACTERISTICS

### Accuracy – Precision

#### Repeatability (Intra-Assay); n = 22

The repeatability was assessed with 2 serum samples under constant parameters (same operator, instrument, day and kit lot).

| Sample | Mean value<br>[U/ml] | CV<br>[%] |
|--------|----------------------|-----------|
| 1      | 19.98                | 2.2       |
| 2      | 3.84                 | 5.0       |

### Reproducibility (Inter-Assay); n = 20

The reproducibility was assessed with 3 serum samples under varying parameters (different operators, instruments, days and kit lots).

| Sample | Mean value<br>[U/ml] | CV<br>[%] |
|--------|----------------------|-----------|
| 1      | 2.98                 | 9.0       |
| 2      | 11.26                | 8.9       |
| 3      | 23.58                | 8.7       |

### Accuracy – Trueness

The trueness states the closeness of the agreement between the result of a measurement and the true value of the measurand. Therefore, DAO-spikes with known concentrations were added to 4 different serum samples. The samples were diluted by the volume of the spike. This was considered when calculating the expected values.

| Sample<br>[U/ml] | Spike<br>[U/ml] | Expected<br>[U/ml] | Obtained<br>[U/ml] | Recovery<br>[%] |
|------------------|-----------------|--------------------|--------------------|-----------------|
| 4.72             | 5.0             | 9.24               | 9.93               | 107.41          |
|                  | 2.5             | 6.98               | 6.89               | 98.64           |
|                  | 1.5             | 6.07               | 5.93               | 97.69           |
| 12.75            | 5.0             | 16.47              | 16.91              | 102.67          |
|                  | 2.5             | 14.61              | 14.21              | 97.24           |
|                  | 1.5             | 13.87              | 13.14              | 94.79           |
| 5.10             | 5.0             | 9.59               | 8.54               | 89.02           |
|                  | 2.5             | 7.34               | 6.59               | 89.75           |
|                  | 1.5             | 6.45               | 6.06               | 93.98           |
| 6.59             | 5.0             | 10.93              | 11.76              | 107.60          |
|                  | 2.5             | 8.76               | 8.89               | 101.52          |
|                  | 1.5             | 7.89               | 7.90               | 100.18          |

### Linearity

The linearity states the ability of a method to provide results proportional to the concentration of analyte in the test sample within a given range. This was assessed according to CLSI guideline

EP06-A with a serial dilution of 5 different serum samples.

For DAO in serum, the method has been demonstrated to be linear from 0.41 to 9.18 U/ml based on the standard curve without considering possibly used sample dilution factors, showing a non-linear behaviour of less than  $\pm 20\%$  in this interval.

| Sample | Dilution | Expected [U/ml] | Obtained [U/ml] | Recovery [%] |
|--------|----------|-----------------|-----------------|--------------|
| 1      | 1:5      | 9.18            | 9.18            | 100.00       |
|        | 1:10     | 4.59            | 4.63            | 100.82       |
|        | 1:20     | 2.29            | 2.29            | 100.04       |
|        | 1:40     | 1.15            | 1.18            | 102.66       |
| 2      | 1:5      | 3.30            | 3.30            | 100.00       |
|        | 1:10     | 1.65            | 1.72            | 104.52       |
|        | 1:20     | 0.82            | 0.88            | 106.58       |
|        | 1:40     | 0.41            | 0.45            | 110.22       |
| 3      | 1:5      | 6.35            | 6.35            | 100.00       |
|        | 1:10     | 3.17            | 3.14            | 98.90        |
|        | 1:20     | 1.59            | 1.88            | 118.59       |
|        | 1:40     | 0.79            | 0.94            | 118.82       |
| 4      | 1:10     | 5.78            | 5.76            | 99.57        |
|        | 1:20     | 2.89            | 3.30            | 113.99       |
|        | 1:40     | 1.45            | 1.79            | 123.58       |
| 5      | 1:10     | 6.92            | 6.82            | 98.57        |
|        | 1:20     | 3.46            | 3.96            | 114.38       |
|        | 1:40     | 1.73            | 2.09            | 120.98       |

### Analytical sensitivity

The following values have been estimated based on the concentrations of the standard without considering possibly used sample dilution factors.

Limit of blank, LoB

0.067 U/ml

|                            |            |
|----------------------------|------------|
| Limit of detection, LoD    | 0.130 U/ml |
| Limit of quantitation, LoQ | 0.195 U/ml |

The evaluation was performed according to the CLSI guideline EP-17-A2. The specified accuracy goal for the LoQ was 20 % CV.

## 12. PRECAUTIONS

- All reagents in the kit package are for *in vitro* diagnostic use only.
- Human materials used in kit components were tested and found to be negative for HIV, Hepatitis B and Hepatitis C. However, for safety reasons, all kit components should be treated as potentially infectious.
- Kit reagents contain sodium azide or ProClin as bactericides. Sodium azide or ProClin are hazardous to health and the environment. Substrates for enzymatic colour reactions may also cause skin and/or respiratory irritation. Any contact with the substances must be avoided. Further safety information can be found in the safety data sheet, which is available from Immundiagnostik AG on request.
- The 10x Wash buffer concentrate (WASHBUF) contains surfactants which may cause severe eye irritation in case of eye contact.  
**Warning:** Causes serious eye irritation. **IF IN EYES:** Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing. If eye irritation persists: get medical Advice/attention.
- The stop solution consists of diluted sulphuric acid, a strong acid. Although diluted, it still must be handled with care. It can cause burns and should be handled with gloves, eye protection, and appropriate protective clothing. Any spill should be wiped up immediately with copious quantities of water. Do not breath vapour and avoid inhalation.

## 13. TECHNICAL HINTS

- Do not interchange different lot numbers of any kit component within the same assay. Furthermore we recommend not assembling wells of different microtiter plates for analysis, even if they are of the same batch.
- Control samples should be analysed with each run.
- Reagents should not be used beyond the expiration date stated on kit label.
- Substrate solution should remain colourless until use.

- To ensure accurate results, proper adhesion of plate sealers during incubation steps is necessary.
- Avoid foaming when mixing reagents.
- Do not mix plugs and caps from different reagents.
- The assay should always be performed according to the enclosed manual.

## 14. GENERAL NOTES ON THE TEST AND TEST PROCEDURE

- This assay was produced and distributed according to the IVD guidelines of 98/79/EC.
- The guidelines for medical laboratories should be followed.
- IDK® is a trademark of Immundiagnostik AG.
- Incubation time, incubation temperature and pipetting volumes of the components are defined by the producer. Any variation of the test procedure, which is not coordinated with the producer, may influence the results of the test. Immundiagnostik AG can therefore not be held responsible for any damage resulting from incorrect use.
- Warranty claims and complaints regarding deficiencies must be logged within 14 days after receipt of the product. The product should be sent to Immundiagnostik AG along with a written complaint.

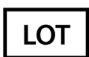













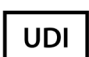

## 15. REFERENCES

1. Sattler J et al. (1988) Food induced histaminosis as an epidemiological problem: plasma histamine elevation and haemodynamic alterations after oral histamine administration and blockade of diamine oxidase (DAO). *Agents and Actions* **23**:361-65.
2. Tufvesson G et al. (1969) Determination of DAO-activity in normal human blood serum. *Scand J Clin Lab Invest* **24**:163-68.
3. Wantke F et al. (1999) The red wine maximization test: drinking histamine rich wine induces a transient increase of plasma diamine oxidase activity in healthy volunteers. *Inflammation Research* **48**:169-70.
4. Wantke F et al. (1994) The red wine provocation test: intolerance to histamine as a model for food intolerance. *Allergy Proceedings* **15**:27-32.



5. Wantke F et al. (1998) Daily variations of serum diamine oxidase and the influence of H1 and H2 blockers: a critical approach to routine diamine oxidase assessment. *Inflammation Research* **47**:396-400.
6. Jarisch R et al. (1999) Role of food allergy and food intolerance in recurrent urticaria. In: Wüthrich B (Hrsg): The Atopy Syndrome in the Third Millennium. *Curr Probl Dermatol*, Basel, Karger, **28**:64-73.
7. Wantke F et al. (1993) Histamine free diet: treatment of choice for histamine induced food intolerance and supporting treatment for chronic headaches *Clin Exp Allergy* **23**: 982-85.
8. Götz M et al. (1996) Histamin-Intoleranz und Diaminoxidasemangel *Allergologie* **9**: 426-30.
9. Jarisch Reinhart, Histamin-Intoleranz. **1. Auflage** (1999), Thieme-Verlag, Stuttgart

**Used symbols:**

|  |  |   |                                    |
|--|--|---|------------------------------------|
|    | Lot number                                 |    | Catalogue number                   |
|    | In Vitro Diagnostic Medical Device         |    | To be used with                    |
|    | Manufacturer                               |    | Contains sufficient for <n> tests  |
|   | Temperature limitation                     |   | Use by                             |
|  | Consult product specification data sheet   |  | Consult instructions for use       |
|  | European Conformity                        |  | Do not re-use                      |
|  | Contains plasma derivatives or human blood |  | Contains material of animal origin |
|  | Unique Device Identification               |  | Irritant                           |

For Reference Use Only

For Reference Use Only

**Immundiagnostik AG**

Stubenwald-Allee 8a  
64625 Bensheim, Germany

Tel.: +49 6251 70190-0

Fax: +49 6251 70190-363

[info@immundiagnostik.com](mailto:info@immundiagnostik.com)

[www.immundiagnostik.com](http://www.immundiagnostik.com)

