

**1. Identification of product and manufacturer**

- 1.1 Product name: MPO, Human, ELISA kit  
Catalog number: **HK324-01, HK324-02**
- 1.2 Intended uses: For research use only. Not for use in or on humans and animals or for diagnostics.
- 1.3 Manufacturer: Hycult Biotech  
Frontstraat 2a  
5405 PB Uden  
The Netherlands  
Tel.: +31 (0)413 251335  
www.hycultbiotech.com  
info@hycultbiotech.com
- 1.4 Emergency phone #:112 in Europe and 911 in America

**2. Hazards identification**

The following hazard Classification and Label elements are attributable to the Stop Solution, which contains 2% Oxalic Acid. The ingredients present in the buffers, standards and detection antibody are in an amount that does not require labeling accordance with EC directives or respective national laws.

**2.1 Classification of the substance or mixture:**

Classification according to Regulation (EC) No 1272/2008

Acute toxicity, Oral (Category 4): **H302**: Harmful if swallowed.

Acute toxicity, Dermal (Category 4), **H312**: Harmful in contact with skin.

Serious eye damage (Category 1), **H318**: Harmful if swallowed or in contact with skin.

**2.2 Label elements**

Labelling according Regulation (EC) No 1272/2008



**Pictogram:**

**Signal Word:** Danger

**Hazard statement(s):**  
**H302**: Harmful if swallowed  
**H312**: Harmful in contact with skin  
**H318**: Causes serious eye damage.

**Precautionary statement(s):** **P305 + P351 + P338; IF IN EYES:** Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing.

**Supplemental Hazard Statements:** none

### **2.3 Other hazards**

The standard is from human origin, which is collected from donors in an FDA-approved collection center. Each unit is tested for the standard FDA-required viral markers, and found negative using FDA-approved methods. While every measure is taken to ensure the safety of our products, we recommend following universal precautions as laid out the following organizations WHO, CDC and OSHA, handle biological materials as if capable of spreading infectious disease.

### **Ecological Information**

This substance/mixture contains no components classified as persistent, bioaccumulative, and toxic (PBT) or very persistent and very bioaccumulative (vPvB) at concentrations of 0.1% or higher.

Additionally, it does not contain components with endocrine-disrupting properties as defined by REACH Article 57(f) at levels of 0.1% or higher.

### **Toxicological Information**

The substance/mixture does not contain components with endocrine-disrupting properties as defined by REACH Article 57(f) at concentrations of 0.1% or higher.

### **3. Information on ingredients**

The following chemical substance can be found in the product, **HK324-01, HK324-02** contains:

| <b>Component</b>                                         | <b>Substance</b>  | <b>Identification number</b>                                     | <b>Concentration<br/>Hazard Limit (%)</b> | <b>Concentration<br/>in %</b> |
|----------------------------------------------------------|-------------------|------------------------------------------------------------------|-------------------------------------------|-------------------------------|
| Buffers,<br>standards<br>and<br>detection<br>antibody    | 2-chloroacetamide | 616-036-00-0 (Index No)<br>201-174-2 (EC No)<br>79-07-2 (CAS No) | ≥ 0.1                                     | 0.02                          |
| Lyophilized<br>standards<br>and<br>detection<br>antibody | 2-chloroacetamide | 616-036-00-0 (Index No)<br>201-174-2 (EC No)<br>79-07-2 (CAS No) | ≥ 0.1                                     | 0.02                          |
| Stop solution                                            | Oxalic acid       | 607-006-00-8(Index No)<br>205-634-3 (EC No)<br>144-62-7 (CAS No) | ≥ 1                                       | 2.0                           |

### **4. First aid measures**

**IF SWALLOWED:** Never give anything by mouth to an unconscious person. Rinse mouth with water.

**IF ON SKIN:** Wash off with soap and plenty of water

**IF IN EYES:** Flush eyes with water as a precaution.

**IN ALL CASES OF DOUBT OR WHEN SYMPTOMS PERSIST, ALWAYS SEEK MEDICAL ATTENTION.**

### **5. Firefighting measures**

Use water spray jet, dry chemicals, carbon dioxide, foam or sand depending on the surrounding equipment and materials. Prevent fire extinguishing water from contaminating surface water or the ground water system.

There are no unusual fire or explosion hazards known for any of the components.

## **6. Accidental release measures**

6.1 Personal precautions: Avoid direct contact with skin, eyes and clothing. Do not inhale vapors or fumes.

6.2 Environmental precautions: Prevent contamination of soil, drains and surface water. Dispose of in accordance with all applicable laws and regulations.

6.3 Cleaning procedure: Soak up with inert absorbent material. After removing the substance, clean the contaminated area and disinfect.

## **7. Handling and storage**

For optimal performance, it is recommended to work according to Good Laboratory Practice (GLP). Avoid prolonged exposure to light. Store in a cool, dry, well-ventilated place in securely closed original packaging. Please refer to labeling and CoA for correct storage and handling of the product.

## **8. Exposure controls / personal protection**

There is no exposure limit value known. The following precautionary measures are recommended as a means of personal protection:

Skin: Wear laboratory protective clothing and closed footwear.

Respiratory protection: None

Eye protection: It is recommended to use safety goggles

Hands: Use protective gloves that are chemically resistant.

Use only with adequate ventilation. Please make sure to wash hands thoroughly after handling chemical products or before eating, smoking and using the restroom.

## **9. Physical and chemical properties**

|                                  |                     |                                      |    |
|----------------------------------|---------------------|--------------------------------------|----|
| <u>Appearance:</u>               | Liquid              |                                      |    |
| <u>Color:</u>                    | Colorless           |                                      |    |
| <u>Odor:</u>                     | Odorless            |                                      |    |
| <u>pH:</u>                       | Neutral             |                                      |    |
| <u>Solubility:</u>               | Soluble in water    |                                      |    |
| <u>Boiling point:</u>            | Not determined (ND) | <u>Decomposition temperature:</u>    | ND |
| <u>Evaporation rate:</u>         | ND                  | <u>Viscosity:</u>                    | ND |
| <u>Vapor pressure:</u>           | ND                  | <u>Vapor density:</u>                | ND |
| <u>Relative density:</u>         | ND                  | <u>Melting point/freezing point:</u> | ND |
| <u>Flash point:</u>              | ND                  | <u>Flammability:</u>                 | ND |
| <u>Autoignition temperature:</u> | ND                  | <u>Particle characteristics:</u>     | ND |

## **10. Stability and reactivity**

Product is stable if handled according to the described storage conditions and before the mentioned expiry date. Use only sterile glass and plastics (polypropylene) suitable for laboratory use.

The products have no known hazardous reactions or decomposition products. Avoid freezing and prolonged exposure to light.

## **11. Toxicological information**

Except for the stop solution, the kit components are in small sizes with a concentration below the acceptable limit for hazardous ingredients. The toxicological risk is minimal. No toxicological experiments have been performed with the kit components.

## **12. Ecological information**

Except for the stop solution, the kit components are in small sizes with a concentration below the acceptable limit for hazardous ingredients. The environmental impact of the kit and all of its components has not been evaluated. There are no significant ecological effects or critical hazards known.

## **13. Disposal considerations**

The generation of waste should be avoided or minimized wherever possible. The product and packaging must be disposed in accordance with all applicable laws and regulations.

## **14. Transport information**

This product is not regulated or classified for transportation of hazardous goods by road, air, rail or sea.

## **15. Regulatory information**

This SDS has been prepared in accordance with the requirements of Regulations (EC) No 1907/2006, (EC) No 1272/2008, (EU) No 2020/878, and (EU) No 2024/2564. For research use only. Not for use in or on humans, animals, or for diagnostic purposes.

## **16. Other information**

Employers must ensure that Safety Data Sheets are readily accessible to employees for all hazardous chemicals in their workplace. All SDS's for Hycult Biotech ELISA kits can be downloaded from the website via the respective product pages.

The information provided is believed to be correct and is intended as a guide to users. Hycult Biotech cannot be held responsible or be liable for the accuracy or completeness of the information contained herein. All materials may present unknown hazards and should be used with caution. Although certain hazards are described herein, we cannot guarantee that these are the only hazards that exist.

Read instructions carefully prior to use. Products may only be handled by technically qualified laboratory personnel. GLP is required.