Myeloperoxidase (human leukocytes)

Haem protein that catalyzes oxidations by H_2O_2 , including MPO-chloride-mediated killing of microbes and tumor cells, inactivation of chemotactic factors, cross-linking of proteins and iodination of proteins.

Ordering Information

Order Online »

| ALX-200-600A-U025 | 25µg |
|-------------------|--------|
| ALX-200-600A-5025 | 5x25µg |

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended. Stored at -20°C

this product is stable for years. Storage of the reconstituted form at 4° C for 2 weeks has no effect on enzyme activity. If solutions must be stored for extended periods of time,

protein concentration should be kept above 1mg/ml and the activity should be

determined prior to use. Freezing solutions of Myeloperoxidase may cause substantial

loss of activity.

Short Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name MPO

Appearance Green solid

CAS 9003-99-0

Formulation Lyophilized from 5mM phosphate buffer, pH 7.0.

Purity ≥98%

Reconstitution Reconstitution of 25 μg with 100 μl distilled water gives

approximately 3.45µM calculated with a molar extinction

coefficient of 91.000 M-1cm-1. per hem at 429nm.

Source Isolated from human leukocytes. Negative for HBsAG-,

HCV- and HIV-antibodies.

UniProt ID P05164

Last modified: May 29, 2024

