## **Granzyme A (human)**

Citations: 3

Ordering Information

Order Online >>

ALX-200-605-C010

10µg

Manuals, SDS & CofA

View Online »

## **Handling & Storage**

**Use/Stability** As indicated on product label or CoA when stored as recommended. Stable for at least

6 months when stored at -80°C. Do not store dilute solutions (1-1000ng/ml) for more

than 24 hours.

**Handling** For maximum product recovery after thawing, centrifuge the vial before opening the cap.

Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.

**Long Term Storage** -80°C

Shipping Dry Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

**Formulation** Liquid. In ~0.15M sodium chloride.

**MW** ~52kDa (fully glycosylated homodimer)

Purity Detail Purified by differential solubilization and cation exchange

chromatography. Dialyzed against 0.15M sodium chloride

and concentrated.

Source Human cytotoxic granules of NK-92 cells.

Specific Activity ~9.2U/µg protein. One unit is defined as the amount of

enzyme that hydrolyzes 1nmole of Z-Lys-SBzl per min. at

25°C in 0.05M TRIS, pH 8.0, containing 0.15M NaCl,

0.01% Triton X-100, 0.3mM DTNB.

UniProt ID P12544

Last modified: May 29, 2024

