Granzyme A (human), (recombinant)

Citations: 1

View Online »

Ordering Information

Order Online »

ALX-201-118-C010

10µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability Stable for at least 6 months after receipt when stored at -80°C. Do not dilute to more

than 0.2mg/ml. 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

Concentration 0.2mg/ml

Formulation Liquid. In 0.05M MES, pH 6.0, containing 10% sucrose

and 0.5M sodium chloride. Contains no preservatives.

MW ~52kDa (homodimer)

Purity ≥98% (SDS-PAGE)

Source Produced in *E. coli*. Active human granzyme A.

Specific Activity Refer to the lot specific CofA for specific activity.

UniProt ID P12544

Last modified: December 17, 2024



eu@enzolifesciences.com