Granzyme B (human) monoclonal antibody (GrB7)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ALX-804-198A-0050

50µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application IHC (PS), WB

Application NotesDetects a band of ~33kDa by Western blot.

Not recommended for IHC (FS).

Clone GrB7

Crossreactivity Does not cross-react with granzyme A.

Formulation Liquid. In PBS containing 0.1% BSA and 0.02% sodium

azide.

Host Mouse

Immunogen Human granzyme B.

lgG2a

Recommendation Dilutions/Conditions Immunohistochemistry (paraffin sections, 1:20)Suggested

dilutions/conditions may not be available for all

applications. Optimal conditions must be determined

individually for each application.

Species Reactivity Human

UniProt ID P10144

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com