Bak polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Citations: 8

View Online »

Ordering Information

Order Online »

ALX-210-002-R050

50µl

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability Stable for at least 1 year after receipt when stored at -80°C.

Long Term Storage -80°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name BCL2L7

Application IHC (PS), WB

Application NotesDetects a band of ~24kDa by Western blot.

Formulation Liquid. In PBS containing 0.05% sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to aa 14-36

(C14GEPALPSASEEQVAQDTEEVFR36) of human Bak.

Recommendation Dilutions/Conditions Western Blot (1-2µl/ml)Suggested dilutions/conditions may

not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity Human, Mouse, Rat

UniProt ID Q16611

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

.

Last modified: May 29, 2024



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com