KDEL monoclonal antibody (10C3) (DyLight™ 488 conjugate)

The tetrapeptide KDEL, located at the carboxy-terminal sequences of luminal proteins, is a retrieval motif essential for the precise sorting of these proteins along the secretory pathway. KDEL proteins perform essential functions in the endoplasmic reticulum (ER) related to protein folding as well as assembly. The localization of chaperones and other soluble proteins to the ER is achieved by their continuous retrieval from post-ER compartments by the KDEL receptor (Erd2p), which is a membrane protein localized in the Golgi apparatus.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

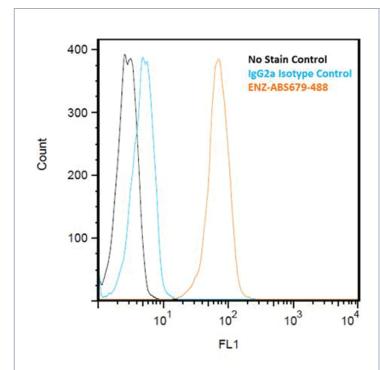
Order Online »

ENZ-ABS679-488-0050

50µg

Manuals, SDS & CofA

View Online »



Flow cytometry analysis of 10⁶ Jurkat cells stained using KDEL monoclonal antibody (10C3), (Prod. No. ENZ-ABS679-488), at a concentration of 10µg/ml.

Handling & Storage

Handling Avoid vigorous centrifugation and vortexing. Do not freeze. Protect from light.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Lys-Asp-Glu-Leu, Lysine-aspartic acid-glutamate-leucine

Application Flow Cytometry

Application NotesDetects three bands of ~94kDa, ~78 kDa and ~55kDa by

Western blot.

Clone 10C3

Formulation Liquid. In PBS containing 0.09% sodium azide.

GenBank ID M14050

Host Mouse

Immunogen Synthetic peptide corresponding to aa 649-654

(S649EKDEL654) of rat Grp78.

Isotype IgG2a

Purity Detail Protein G affinity purified.

Recommendation Dilutions/Conditions Flow Cytometry (1:20)Suggested dilutions/conditions may

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

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Avian, Insect, Mammalian, Plant, Yeast

Specificity Recognizes proteins containing the KDEL sequence.

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-

eu@enzolifesciences.com