

Data Sheet

Anti-ERK1/2

Mouse Monoclonal Antibody
Anti-ERK1/2 (E31R) Signal Transduction
Catalog #: 25006

Specificity:

Recognizes human ERK1 and ERK2 proteins

Immunogen:

Full length recombinant human ERK1 protein

Applications:

WB

Image:

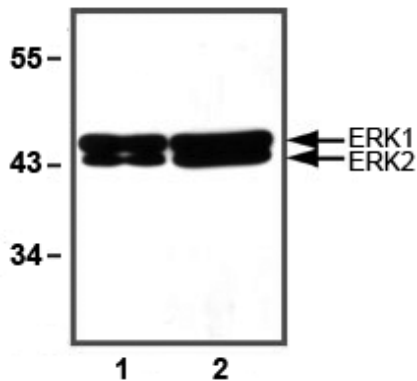


Fig. 1:1000 (1µg/mL) Ab dilution used in WB of 10µg (1) and 30µg (2) HeLa cell lysates

Clonality | Host Species | Isotype:

E31R | Mouse | IgG1

Formulation:

1mg/mL in 10mM PBS, pH 7.2, with 0.05% NaN₃

Cross Reactivity:

WB: ERK1 and ERK2 from human cells. ERK1 and ERK2 from other species may also be detectable

Working Conditions:

WB (with ECL): 1:1000 dilution (incubate for one hour at room temperature)
For best results with other assays, please determine optimal working dilution by titration test

Purification:

Protein A affinity chromatography from mouse ascites fluid

Storage and Stability:

Store at -20°C for up to 2 years. Centrifuge after first thaw to maximize product recovery. Aliquot to avoid repeated freeze-thaw cycles. Store aliquots at 4°C for 1-2 weeks

OUR PRODUCTS ARE FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com