

Data Sheet

Anti-DYKDDDDK (FLAG®) Epitope Tag) Antibody, Monoclonal

Mouse Monoclonal Antibody
Anti-FLAG Tag (FG4R)
Catalog #: 25003 Lot#: 220315

Specificity:

Recognizes the N-terminal, C-terminal or internal FLAG-tagged fusion proteins

Immunogen:

DYKDDDDK (FLAG) synthetic peptide conjugated to KLH

Applications:

Dot Blot, ELISA, IP, IS, WB

image:

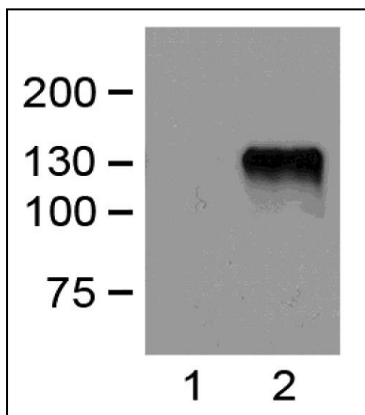


Fig. 1:1000 (1 µg/mL) Ab dilution probed against HEK293 cells transfected with FLAG-tagged protein vector; untransfected (1) and transfected (2).

Clonality | Host Species | Isotype:

FG4R | Mouse | IgG1

Formulation:

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

Sensitivity:

WB (with ECL): 1-5 ng of FLAG-tagged fusion proteins

Working Conditions:

WB (with ECL): 1:1000-3000 dilution (incubate for one hour at room temperature)

IP: 1:50-200 dilution

IS: 1:500-2000 dilution

For best results with other assays (e.g.: Dot, ELISA, etc), please determine optimal working dilution by titration test

Purification:

Protein A affinity chromatography from mouse ascites fluid

Storage:

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 **1.858.202.1401** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com