

FBXO6 Monoclonal Antibody [3F10]

[Catalog No. A-8127-100]

Description

Mouse monoclonal antibody raised against a full length recombinant FBXO6, clone 3F10.

Size

100 µg

Purification

Protein A purified

Immunogen

FBXO6 (AAH20880, 1 a.a. ~ 294 a.a) full length recombinant protein with GST tag. MW of the GST tag alone is 26 KDa.

Cross Reactivity

Human

Isotype

IgG1 Kappa

Formulation

In 1x PBS, pH 7.2

Storage

Aliquot and store at -20°C or lower. Avoid repeated freezing and thawing. Multiple freeze/thaw cycles may result in decreased performance. Non-hazardous. No MSDS necessary.

Handling Recommendations

For maximum recovery of the products, centrifuge the vial prior to opening the cap.

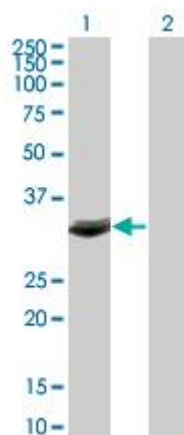
Applications & Suggested Dilutions*

Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

**The end user is responsible for determining optimal working dilutions.*

Data



Western Blot analysis of FBXO6 expression in transfected 293T cell line by FBXO6 Monoclonal Antibody, clone 3F10.

Lane 1: FBXO6 transfected lysate (33.933 KDa).

Lane 2: Non-transfected lysate.

This product is for research purposes only. Not intended for use in diagnostic procedures.