
FBXO24 Monoclonal Antibody [7F12]

[Catalog No. A-8123-100]

Description

Mouse monoclonal antibody raised against a partial recombinant FBXO24, clone 7F12.

Size

100 µg

Purification

Protein A purified

Immunogen

FBXO24 (NP_277041, 1 a.a. ~ 101 a.a) partial 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*

ELISA: 0.25-1 µg/ml

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

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