

FBXO9 Monoclonal Antibody [1G12]

[Catalog No. A-8125-100]

Description

Mouse monoclonal antibody raised against a partial recombinant FBXO9, clone 1G12.

Size

100 µg

Purification

Protein A purified

Immunogen

FBXO9 (NP_036479, 338 a.a. ~ 448 a.a) partial recombinant protein with GST tag. MW of the GST tag alone is 26 KDa.

Cross Reactivity

Human

Isotype

IgG2a 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.

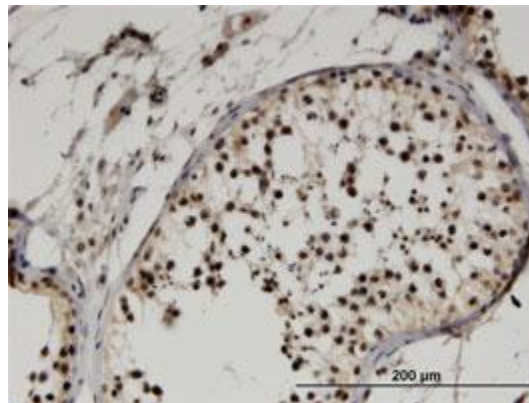
Applications & Suggested Dilutions*

ELISA: 0.25-1 µg/ml

Immunohistochemistry (IHC): 2.5 µg/ml

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

Data



Immunoperoxidase of monoclonal antibody to FBXO9 on formalin-fixed paraffin-embedded human testis. [Antibody Concentration: 3 µg/ml]

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