

CTGF Monoclonal Antibody [2C11]

[Catalog No. A-6216-100]

Description

Mouse monoclonal antibody raised against a full-length recombinant CTGF, clone 2C11.

Size

100 µg

Purification

Protein A purified

Immunogen

CTGF (NP_001892, 31 a.a. ~ 251 a.a) full-length recombinant protein with GST tag. MW of the GST tag alone is 26 KDa.

Cross Reactivity

Human

Isotype

IgG2b 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.

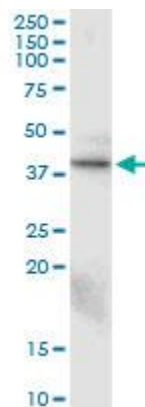
Applications & Suggested Dilutions*

Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

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

Data



CTGF Monoclonal Antibody, clone 2C11. Western Blot analysis of CTGF expression in human kidney.

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