
MBD2 Protein

(Catalog No. E42001)

Description

Human MBD2 (AAH24021.1, 150 a.a. - 220 a.a.) recombinant protein with GST-tag at N-terminal

Molecular Weight (kDa)

34.8 kDa

Formulation

50 mM Tris-HCl, 500 mM sodium chloride, 5% glycerol, and 5 mM β -mercaptoethanol in the elution buffer.

Source

In vitro *E. coli* expression system

Size

10 μ g

Storage Temp & Conditions

Store at -80°C . Aliquot to avoid repeated freezing and thawing. Best use within three months from the date of receipt of this protein.

Purification Notes

>95% in purity

Application Notes

Enzyme-linked Immunoabsorbent Assay; Western Blot (Recombinant protein); Antibody Production; Protein Array.

Ordering Information:

Products	Size	Cat. No.
MBD2 Protein	10 μ g	E42001 -1

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