

MDM2 Protein

Recombinant human protein expressed in E. coli cells

Catalog # M45-31G

Lot # H294-4

Product Description

Recombinant human MDM2 (202-end) was expressed in E. coli cells using an N-terminal GST tag. The gene accession number is [NM_002392](#).

Gene Aliases

HDMX, hdm2, MGC71221

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

Scientific Background

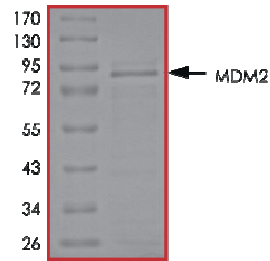
MDM2 is a nuclear phosphoprotein that binds and inhibits transactivation by p53, as part of an autoregulatory negative feedback loop (1). Overexpression of the MDM2 gene product can lead to excessive inactivation of p53 thereby diminishing its tumor suppressor function. The inactivation of p53 is mediated by the E3 ubiquitin ligase activity of MDM2 which targets p53 for proteasomal degradation. MDM2 also affects the cell cycle, apoptosis, and tumorigenesis through interactions with other proteins, including retinoblastoma 1 and ribosomal protein L5 (2). Amplification of MDM2 is frequently observed in human sarcomas and this is consistent with the hypothesis that MDM2 binds to p53 which then leads to escape from p53-regulated growth control.

References

1. Engel, T. et al: Elevated p53 and lower MDM2 expression in hippocampus from patients with intractable temporal lobe epilepsy. *Epilepsy Res.* 2007;77(2-3):151-6
2. Liu, G. et al: Genetic polymorphisms of MDM2, cumulative cigarette smoking and nonsmall cell lung cancer risk. *Int J Cancer.* 2008;122(4):915-8

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

Purity



The purity was determined to be **>85%** by densitometry. Approx. MW **84kDa**.

MDM2 Protein

Recombinant human protein expressed in E. coli cells

Catalog Number	M45-31G
Specific Lot Number	H294-4
Purity	>85%
Concentration	0.2 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.