

MDM4 (1-137) Protein

Recombinant protein expressed in E.coli cells

Catalog # M47-31BH

Lot # A1437-3

Product Description

Recombinant human MDM4 (1-137) was expressed in E. coli cells using an N-terminal His tag. The gene accession number is [NM_002393](#).

Gene Aliases

DKFZp781B1423; HDMX; MDMX; MGC132766; MRP1

Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM DTT, 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.

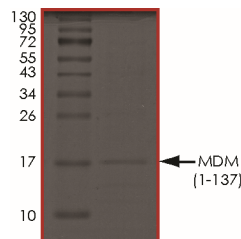
Scientific Background

MDM4 is a nuclear protein that contains a p53 binding domain at the N-terminus and a RING finger domain at the C-terminus. MDM4 shows structural similarity to p53-binding protein MDM2 and both proteins bind the p53 tumor suppressor protein and inhibit its activity. However, unlike MDM2, MDM4 does not cause nuclear export or degradation of p53. Instead, MDM4 inhibits p53 activity by binding to the transcriptional activation domain of p53. MDM4 is overexpressed in a variety of human cancers (1). Expression level of MDM4 is significantly higher in chronic lymphocytic leukemia. MDM4 is a specific chemotherapeutic target for treating retinoblastoma (2).

References

1. Parant, J.et.al: Rescue of embryonic lethality in Mdm4-null mice by loss of Trp53 suggests a nonoverlapping pathway with MDM2 to regulate p53. Nature Genet. 29: 92-95, 2001.
2. Laurie, N. A. et.al: Inactivation of the p53 pathway in retinoblastoma. Nature 444: 61-66, 2006.

Purity



The purity of MDM4 (1-137) was determined to be **>75%** by densitometry.
Approx. MW **17.3 kDa**.

MDM4 (1-137) Protein

Recombinant protein expressed in E. coli cells

Catalog #	M47-31BH
Lot #	A1437-3
Purity	>75%
Concentration	0.2µg/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.

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

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.