

SMAD5 Protein

Full-length recombinant protein expressed in E. coli cells

Catalog # S14-30G

Lot # E203-3

Product Description

Recombinant full-length human SMAD5 was expressed in E. coli cells using an N-terminal GST tag. The gene accession number is [NM_005903](#).

Gene Aliases

Dwfc, JV5-1, MADH5, DKFZp781C1895, DKFZp781O1323

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.

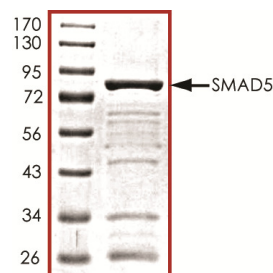
Scientific Background

SMAD5 is a member of the SMAD family and mediates signaling by the transforming growth factor-beta (TGFβ) superfamily and related ligands (1). SMAD5 plays a critical role in the signaling pathway by which TGFβ inhibits the proliferation of human hematopoietic progenitor cells. SMAD5 is up-regulated in gastric epithelial cells during the infection of the pathogen Helicobacter pylori and it mediates apoptosis of gastric epithelial cells induced by H. pylori infection (2). In mature human B cells, bone morphogenetic protein 6 (BMP-6) inhibits cell growth and rapidly induces phosphorylation of SMADs 5 and 8.

References

- Heldin, C.H. et al: TGF-beta signalling from cell membrane to nucleus through SMAD proteins". Nature, 1997, **390** (6659): 465-71.
- Nagasko, T. et al: Up-regulated Smad5 mediates apoptosis of gastric epithelial cells induced by Helicobacter pylori infection. J Biol Chem. 2003 Feb 14;278(7):4821-5.

Purity



The purity of SMAD5 was determined to be **>80%** by densitometry. Approx. MW **82kDa**.

SMAD5 Protein

Full-length recombinant protein expressed in E. coli cells

Catalog Number **S14-30G**

Specific Lot Number **E203-3**

Purity	>80%
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.

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.