

TGFβR1, Unactive

Recombinant protein expressed in Sf9 cells

Catalog # T07-35G

Lot # G071-2

Product Description

Recombinant human TGFβR1 (80-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC071181](#).

Gene Aliases

ST AAT5, ALK5, SKR4, ALK-5, TGFR-1, ACVRLK4

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

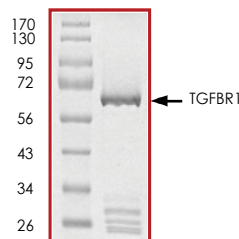
TGFβR1 is a transmembrane serine/threonine protein kinase and a member of the TGFβ receptor subfamily (1). TGFβ regulates cell cycle progression by a unique signaling mechanism that involves its binding to TGFβR2 and activation of TGFβR1. TGFβR1 may be inactivated in many of the cases of human tumor cells refractory to TGFβ-mediated cell cycle arrest. Heterozygous mutations in TGFβR1 and TGFβR2 have been reported in Loeys-Dietz aortic aneurysm syndrome (LDS) and also dominant TGFβR2 mutations have been identified in Marfan syndrome type 2 (MFS2) and familial thoracic aortic aneurysms and dissections (TAAD). Mutations of TGFβR1 and TGFβR2 are associated with atherosclerosis and several human cancers (2).

References

1. Mátyás, G. et al: Identification and in silico analyses of novel TGFβR1 and TGFβR2 mutations in Marfan syndrome-related disorders. Hum Mutat. 2006;27(8):760-9.
2. Suarez, B.K. et al: TGFβR1*6A is not associated with prostate cancer in men of European ancestry. Prostate Cancer Prostatic Dis. 2005;8(1):50-3.

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 of TGFβR1 was determined to be **>90%** by densitometry. Approx. MW **65kDa**.

TGFβR1, Unactive

Recombinant protein expressed in Sf9 cells

Catalog Number	T07-35G
Specific Lot Number	G071-2

Purity	>90%
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.