

TGFβR1 Peptide

Synthetic peptide substrate derived from human SMAD3

Catalog # T36-58

Lot # 1087-2

Product Description

The TGFβR1 peptide sequence (KKKVLQMGSPSIRC-S(pS)VS) is derived from human SMAD3 (215-230) and is suitable for use as the substrate for TGFβR1 superfamily, including ACVRs (ALK1, ALK2, ALK4 and ALK7) and BMPRs (ALK3 and ALK6).

Molecular Weight

The theoretical molecular weight is 2116.45

Purity

The purity was determined to be 98.1% by HPLC analysis.

Formulation

1mg of peptide supplied as a lyophilized powder.

Storage and Stability

Store product at -20°C. For optimal storage, aliquot diluted product into smaller quantities and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

Reconstitution Protocol

Dilute peptide in 20mM Tris-HCl, pH 7.5 solution to a final concentration of 1mg/ml.

Related Kinases

TGFβR1 Peptide can be utilized as a substrate for the following active protein kinases.

Product Name	Catalog Number
ALK1	A09-11G
ALK2	A06-11G
ALK4	A07-11G

TGFβR1 Peptide

Synthetic peptide substrate derived from human SMAD3

Catalog #	T36-58
Quantity	1 mg
Lot #	1087-2
Purity	98.1%
Format	1 mg lyophilized powder
Stability	1yr at -20°C from date of shipment
Storage & Shipping	Store product at -20°C. For optimal storage, aliquot diluted product into smaller quantities and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped at ambient temperature.