

TGF β R1 Peptide

Synthetic peptide substrate derived from human SMAD3

Catalog # T36-58

Lot # E105-4

Product Description

The TGF β R1 peptide sequence (KKKVLQMGSPSIRC-S(pS)VS) is derived from human SMAD3 (215-230) and is suitable for use as the substrate for TGF β R1 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 95.79% 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

TGF β R1 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

TGF β R1 Peptide

Synthetic peptide substrate derived from human SMAD3

Catalog Number	T36-58
Quantity	1 mg
Specific Lot Number	E105-4
Purity	95.79%
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.