

Product Data Sheet

Product Name: TIF2 (740-753), Transcriptional Intermediary Factor 2 (740-753)

Catalog Number: AS-61992 (1 mg) Lot Number: See label on vial

Sequence: H-Lys-Glu-Asn-Ala-Leu-Leu-Arg-Tyr-Leu-Leu-Asp-Lys-Asp-Asp-OH

(3-letter code)

KENALLRYLLDKDD (1-letter code)

Molecular Weight: 1706.9 % Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Using H₂O, reconstitute by adding 100 μl to 1 mg TIF2 peptide.

Storage: TIF2 peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at –20°C or lower. Reconstituted peptide can be aliquoted and stored at –20°C or lower.

Description: This peptide is a nuclear receptor (NR) box B3 region of the p160 co-activator Transcriptional Intermediary Factor 2 (TIF2) peptide, a LXXLL motif. The activation function 2/ligand-dependent interaction between nuclear receptors and their co-regulators is mediated by a short consensus motif nuclear receptor box. Wärnmark, A., et al. *J. Biol. Chem.* **277**, 21862 (2002).

Additional Information: Listed below are relevant information that may provide a guideline on how to use this product. End users will have to adapt to their own specific applications.

Both TIF2 and NCoR peptides were obtained from AnaSpec Inc. SPR binding surfaces were prepared by immobilizing approximately 300–400 RU of biotinylated coactivator (TIF2) or corepressor (NCoR-ID2) peptides on NeutrAvidin charged CM4 chips. Solution competition experiments were performed by incubating a range of TIF2 or NCoR-ID2 peptide concentrations with 200 nM GR-LBD(F602S) bound with Dexamethasone or RU-486. These reaction mixtures were then injected over immobilized TIF2 and NCoR-ID2 surfaces at a flow rate of 30 μL/min until equilibrium was reached-Kroe, RR. et al. *Biophys. Chem.* 128, 156 (2007).

Published Citations:

Kroe, RR. et al. Biophys. Chem. 128, 156 (2007).

For Research Use Only