



## Product Data Sheet

---

<b>Product Name:</b>	Insulin Receptor (1142-1153), pTyr1150	
<b>Catalog Number:</b>	AS-20274 (1 mg)	Lot Number: See label on vial
<b>Sequence:</b>	H-Thr-Arg-Asp-Ile-Tyr-Glu-Thr-Asp-pTyr-Tyr-Arg-Lys-OH (3-letter code) TRDIYETD-pY-YRK (1-letter code)	
<b>Molecular Weight:</b>	1702.7	
<b>Peptide Purity:</b>	>95%	
<b>Appearance:</b>	Lyophilized white powder	

**Peptide Reconstitution:** Insulin Receptor (1142-1153) peptide is freely soluble in water.

**Storage:** Insulin Receptor (1142-1153) peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at  $-20^{\circ}\text{C}$  or lower. Reconstituted peptide can be aliquoted and stored at  $-20^{\circ}\text{C}$  or lower.

**Description:** Peptides are used as insulin receptor tyrosine kinase substrates. Ref: Keane, N. et al. *Eur. J. Biochem.* **226**, 525 (1994); Stadtmauer, L. and O. Rosen, *J. Biol. Chem.* **261**, 3402 (1986).

### Related Products:

<b>Name</b>	<b>Cat #</b>	<b>Size</b>
Insulin Receptor (1142-1153), pTyr1146 TRDI-pY-ETDYRK	AS-20292	1 mg

*For Research Use Only*