



## Product Data Sheet

---

<b>Product Name:</b>	Prosaptide TX14(A)	
<b>Catalog Number:</b>	AS-60248-1 (1 mg) AS-60248-5 (5 mg)	Lot Number: See label on vial
<b>Sequence:</b>	H-Thr-D-Ala-Leu-Ile-Asp-Asn-Asn-Ala-Thr-Glu-Glu-Ile-Leu-Tyr-OH (3-letter code) TaLIDNNATEEILY (1-letter code)	
<b>Molecular Weight:</b>	1579.7	
<b>Peptide Purity:</b>	>95%	
<b>Appearance:</b>	Lyophilized white powder	

**Peptide Reconstitution:** Prosaptide TX14(A) peptide is freely soluble in 1% NH<sub>4</sub>OH.

**Storage:** Prosaptide TX14(A) 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 14-mer prosaptide sequence is derived from the active neurotrophic region in the amino-terminal portion of the saposin C domain. Synthetic peptides derived from this region are biologically active and are named "prosaptides." Prosaposin and prosaptides are active on a variety of neuronal cells, stimulating sulfatide synthesis and increasing sulfatide concentration in Schwann cells and oligodendrocytes. This indicates that prosaposin and prosaptides are trophic factors for myelin formation. Ref: Campana, M. et al. *FASEB J.* **12**, 307 (1998); Hiraiwa, M. et al. *Proc. Natl. Acad. Sci. USA* **94**, 4778 (1997); Taylor, E. et al. *Pharmacol.* **295**, 190 (2000); Calcutt, NA. et al. *Anesthesiology* **93**, 1271 (2000).

**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.

TX14(A), a prosaptide (TX14(A)=TXLIDNNATEEILY, where X equals D-alanine) derived from the neurotrophic region of saposin C, was synthesized commercially to 98% purity (AnaSpec, San Jose, Calif.). iSC cells (approximately 2.0x10<sup>7</sup>) were incubated in DMEM/F12 without serum 18 h before stimulation with TX14(A) for 5 min at 37°C-[Campana, W. M. et al. FASEB J. 12, 307 \(1998\).](#)

### Published Citations:

Hiraiwa, M. et al. *PNAS.* **94**, 4778 (1997).

Campana, W. M. et al. *FASEB J.* **12**, 307 (1998).

Calcutt, NA. et al. *Anesthesiology* **93**, 1271 (2000).. *For Research Use Only*