



## Product Data Sheet

---

<b>Product Name:</b>	Thrombin Receptor Activator for Peptide 6 (TRAP-6)	
<b>Catalog Number:</b>	AS-24190 (1 mg) AS-24191 (5 mg) AS-24190-25 (25 mg)	<b>Lot Number:</b> See label on vial
<b>Sequence:</b>	H-Ser-Phe-Leu-Leu-Arg-Asn-OH (3-letter code) SFLLRN (1-letter code)	
<b>Molecular Weight:</b>	748.9	
<b>Peptide Purity:</b>	>95%	
<b>Appearance:</b>	Lyophilized white powder	

**Peptide Reconstitution:** Initially try to dissolve a small amount of Thrombin Receptor Activator peptide in water. If peptide is insoluble in water, use DMSO.

**Storage:** Thrombin Receptor Activator 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.

**References:** Scarborough, RM. et al. *J. Biol. Chem.* **267**, 13146 (1992); Ahmad, S. and PN. Walsh, *Biochem.* **44**, 13858 (2005); Buegler, M. et al. *J. Thromb. Thrombolysis* **19**, 115 (2005).

### Published Citations:

- Birukova, AA. et al. *FASEB J.* **18**, 1879 (2004).
- Kawkitinarong, K. et al. *Am. J. Respir. Cell Mol. Biol.* **31**, 517 (2004).
- Ludeman, MJ. et al. *J. Biol. Chem.* **279**, 18592 (2004).
- Jenkins, R. et al. *J. Clin. Invest.* **116**, 1606 (2006).

### Related Products:

<b>Name</b>	<b>Cat #</b>	<b>Size</b>
Scrambled TRAP Fragment SFLLRN-NH2	AS-60197-1	1 mg
Thrombin Receptor Agonist, amide SFLLR-NH2	AS-60680	1 mg

For Research Use Only

