

Product Data Sheet

Pannexin-1 (Panx1), Mimetic Blocking Peptide	
AS-61911 (1 mg)	Lot Number: See label on vial
H-Trp-Arg-GIn-Ala-Ala-Phe-Val-Asp-Ser-Tyr-OH (3-letter code) WRQAAFVDSY (1-letter code)	
1242.4	
>95%	
Lyophilized white powder	
	AS-61911 (1 mg) H-Trp-Arg-Gln-Ala-Ala-Phe-Val-Asp WRQAAFVDSY (1-letter code) 1242.4 >95%

Peptide Reconstitution: Pannexin-1 peptide is freely soluble in water.

Storage: Pannexin-1 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 is a Pannexin-1 (Panx1) mimetic blocking peptide. Pannexin-1 is a recently identified membrane protein that can form gap junction-like connections allowing intercellular passage of dyes when overexpressed in two adjacent oocytes or mammalian epithelial cell lines. Blockade of pannexin-1 in macrophage endogenously expressing the ATP-gated P2X7 receptor (P2X7R) blocks the initial dye uptake, but not the ionic current, and also blocks processing and release of interleukin-1 β (IL-1 β) in response to P2X7R activation. Ref: Pelegrin, P. et al. *EMBO. J.* **25**, 5071 (2006); Pelegrin, P. et al.t *J. Biol. Chem.* **282**, 2386 (2007).

For Research Use Only