

# PRODUCT INFORMATION

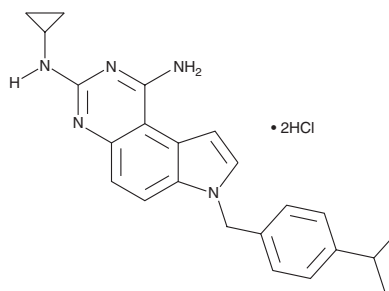


## SCH 79797 (hydrochloride)

Item No. 14633

**CAS Registry No.:** 1216720-69-2  
**Formal Name:** N<sup>3</sup>-cyclopropyl-7-[[4-(1-methylethyl)phenyl]methyl]-7H-pyrrolo[3,2-f]quinazoline-1,3-diamine, monohydrochloride

**MF:** C<sub>23</sub>H<sub>25</sub>N<sub>5</sub> • 2HCl  
**FW:** 444.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 214, 246, 340 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

SCH 79797 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the SCH 79797 (hydrochloride) in the solvent of choice. SCH 79797 (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of SCH 79797 (hydrochloride) in these solvents is approximately 1, 11, and 5 mg/ml, respectively.

SCH 79797 (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SCH 79797 (hydrochloride) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. SCH 79797 (hydrochloride) has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

SCH 79797 is a non-peptide antagonist of proteinase-activated receptor 1 (PAR1).<sup>1</sup> It blocks binding of the high affinity thrombin receptor-activating peptide (haTRAP) (IC<sub>50</sub> = 70 nM) as well as platelet aggregation induced by either haTRAP or α-thrombin (IC<sub>50</sub> = 0.3 and 3 μM, respectively).<sup>2</sup> SCH 79797 has no activity against PAR2, PAR4, or other receptors involved in platelet activation.<sup>2</sup> It also blocks PAR1 activation on vascular smooth muscle cells and endothelial cells.<sup>2,3</sup>

### References

1. Ahn, H.-S., Arik, L., Boykow, G., *et al.* Structure-activity relationships of pyrroloquinazolines as thrombin receptor antagonists. *Bioorg. Med. Chem. Lett.* **9(14)**, 2073-2078 (1999).
2. Ahn, H.-S., Foster, C., Boykow, G., *et al.* Inhibition of cellular action of thrombin by N<sup>3</sup>-cyclopropyl-7-[[4-(1-methylethyl)phenyl]methyl]-7H-pyrrolo[3,2-f]quinazoline-1,3-diamine (SCH 79797), a nonpeptide thrombin receptor antagonist. *Biochem. Pharmacol.* **60(10)**, 1425-1434 (2000).
3. Lidington, E.A., Steinberg, R., Kinderlerer, A.R., *et al.* A role for proteinase-activated receptor 2 and PKC-ε in thrombin-mediated induction of decay-accelerating factor on human endothelial cells. *Am. J. Physiol. Cell Physiol.* **289(6)**, 1437-1447 (2005).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/01/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM