

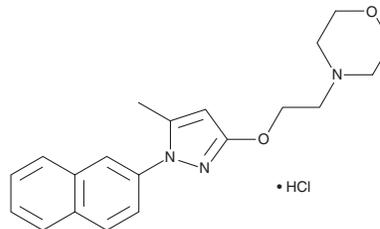
# PRODUCT INFORMATION



## S1RA (hydrochloride)

Item No. 32729

**CAS Registry No.:** 1265917-14-3  
**Formal Name:** 4-[2-[[5-methyl-1-(2-naphthalenyl)-1H-pyrazol-3-yl]oxy]ethyl]-morpholine, monohydrochloride  
**Synonyms:** E-52862, P 027  
**MF:** C<sub>20</sub>H<sub>23</sub>N<sub>3</sub>O<sub>2</sub> • HCl  
**FW:** 373.9  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 217, 258 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

S1RA (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the S1RA (hydrochloride) in the solvent of choice, which should be purged with an inert gas. S1RA (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of S1RA (hydrochloride) in these solvents is approximately 1 and 10 mg/ml, respectively.

S1RA (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the crystalline solid should first be dissolved in DMF and then diluted with the aqueous buffer of choice. S1RA (hydrochloride) has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

S1RA is a sigma-1 ( $\sigma_1$ ) receptor antagonist ( $K_i = 17$  nM).<sup>1</sup> It is selective for  $\sigma_1$  receptors over  $\sigma_2$  receptors ( $K_i = 9,300$  nM), as well as a panel of 170 additional receptors at 1  $\mu$ M. S1RA reduces capsaicin-induced mechanical allodynia, formalin-induced licking and biting behavior, and partial sciatic nerve ligation-induced thermal hyperalgesia in mice ( $ED_{50}$ s = 26.3, 43.7, and 18.8 mg/kg, respectively). It also enhances fentanyl- or loperamide-induced analgesia without affecting gastrointestinal transit in mice.<sup>2</sup>

### References

1. Romero, L., Zamanillo, D., Nadal, X., *et al.* Pharmacological properties of S1RA, a new sigma-1 receptor antagonist that inhibits neuropathic pain and activity-induced spinal sensitization. *Br. J. Pharmacol.* **166**(8), 2289-2306 (2012).
2. Sánchez-Fernández, C., Montilla-García, Á., González-Cano, R., *et al.* Modulation of peripheral  $\mu$ -opioid analgesia by  $\sigma_1$  receptors. *J. Pharmacol. Exp. Ther.* **348**(1), 32-45 (2014).

#### 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, 11/09/2022

#### 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