

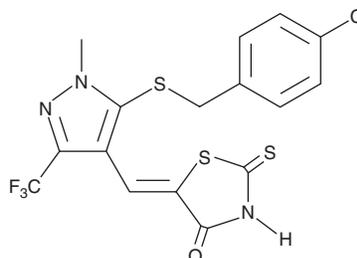
PRODUCT INFORMATION



ADAMTS-5 Inhibitor

Item No. 17188

CAS Registry No.: 929634-33-3
Formal Name: 5-[[5-[[[4-chlorophenyl)methyl]thio]-1-methyl-3-(trifluoromethyl)-1H-pyrazol-4-yl]methylene]-2-thioxo-4-thiazolidinone
MF: C₁₆H₁₁ClF₃N₃OS₃
FW: 449.9
Purity: ≥95%
UV/Vis.: λ_{max}: 225, 280, 363 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

ADAMTS-5 inhibitor is supplied as a crystalline solid. A stock solution may be made by dissolving the ADAMTS-5 inhibitor in the solvent of choice, which should be purged with an inert gas. ADAMTS-5 inhibitor is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ADAMTS-5 inhibitor in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

ADAMTS-5 inhibitor is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ADAMTS-5 inhibitor should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. ADAMTS-5 inhibitor has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

ADAMTS-5 Inhibitor is an inhibitor of ADAMTS-5 (aggrecanase-2; IC₅₀ = 1.1 μM) with 40-fold selectivity over ADAMTS-4 (aggrecanase-1).^{1,2}

References

1. Bursavich, M.G., Gilbert, A.M., Lombardi, S., *et al.* Synthesis and evaluation of aryl thioxothiazolidinone inhibitors of ADAMTS-5 (Aggrecanase-2). *Bioorg. Med. Chem. Lett.* **17(5)**, 1185-1188 (2007).
2. Gilbert, A.M., Bursavich, M.G., Lombardi, S., *et al.* 5-((1H-pyrazol-4-yl)methylene)-2-thioxothiazolidin-4-one inhibitors of ADAMTS-5. *Bioorg. Med. Chem. Lett.* **17(5)**, 1189-1192 (2007).

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, 12/06/2022

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