

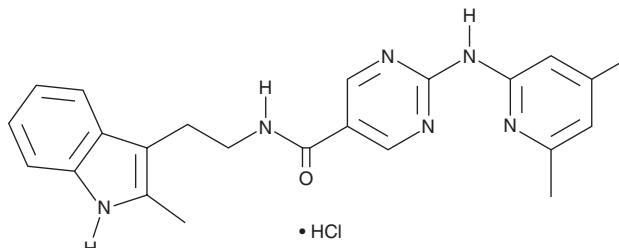
PRODUCT INFORMATION



TG11-77 (hydrochloride)

Item No. 30188

CAS Registry No.: 2550393-38-7
Formal Name: 2-((4,6-dimethylpyridin-2-yl)amino)-N-(2-(2-methyl-1H-indol-3-yl)ethyl)pyrimidine-5-carboxamide, monohydrochloride
MF: $C_{23}H_{24}N_6O \cdot HCl$
FW: 436.9
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 225, 261, 324 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 2 years



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

Laboratory Procedures

TG11-77 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the TG11-77 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. TG11-77 (hydrochloride) is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

TG11-77 (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, TG11-77 (hydrochloride) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. TG11-77 (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

TG11-77 is a brain-permeable antagonist of the prostaglandin E_2 (PGE_2) receptor subtype EP_2 ($K_B = 9.7$ nM).¹ It is selective for EP_2 over DP_1 , EP_4 , and IP receptors ($K_{Bs} = 7,320$, $5,300$, and $>10,000$ nM, respectively). TG11-77 (0.3 and 1 μM) inhibits the expression of genes encoding COX-2, IL-1 β , and IL-6 induced by LPS and the EP_2 receptor agonist ONO-AE1-259-01 in mouse BV-2 microglia expressing human EP_2 receptors.

Reference

1. Amaradhi, R., Banik, A., Mohammed, S., *et al.* Potent, selective, water soluble, brain-permeable EP_2 receptor antagonist for use in central nervous system disease models. *J. Med. Chem.* **63**(3), 1032-1050 (2020).

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, 09/02/2021

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