

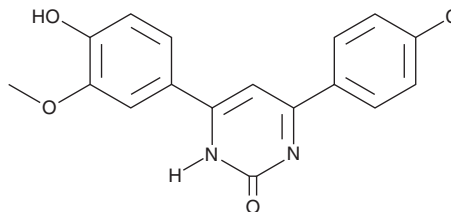
PRODUCT INFORMATION



LIT-927

Item No. 28436

CAS Registry No.: 2172879-52-4
Formal Name: 4-(4-chlorophenyl)-6-(4-hydroxy-3-methoxyphenyl)-2(1H)-pyrimidinone
MF: C₁₇H₁₃ClN₂O₃
FW: 328.8
Purity: ≥98%
UV/Vis.: λ_{max}: 263, 365 nm
Supplied as: A 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

LIT-927 is supplied as a solid. A stock solution may be made by dissolving the LIT-927 in the solvent of choice, which should be purged with an inert gas. LIT-927 is soluble in the organic solvent dimethyl formamide at a concentration of approximately 1 mg/ml.

Description

LIT-927 is an orally bioavailable neutraligand that binds to chemokine (C-X-C motif) ligand 12 (CXCL12; K_i = 267 nM in a FRET-based binding assay).¹ It is selective for inhibiting calcium accumulation induced by CXCL12 over that induced by chemokine (C-C motif) ligand 17 (CCL17), CCL22, CCL5, and CCL2 at 10 μM. It inhibits pulmonary artery smooth muscle cell and pericyte migration *in vitro*.² LIT-927 (100 mg/kg per day) inhibits increases in serum levels of CXCL12, as well as reduces total pulmonary vascular resistance and remodeling of pulmonary arterioles, in a rat model of pulmonary hypertension induced by monocrotaline (Item No. 16666). It inhibits allergen-induced eosinophil infiltration into the lung by 62% in a mouse model of allergic airway hyper eosinophilia when administered at an oral dose of 1,400 μmol/kg.

References

1. Regenass, P., Abboud, D., Daubeuf, F., *et al.* Discovery of a locally and orally active CXCL12 neutraligand (LIT-927) with anti-inflammatory effect in a murine model of allergic airway hyper eosinophilia. *J. Med. Chem.* **61**(17), 7671-7686 (2018).
2. Bordenave, J., Thuillet, R., Tu, L., *et al.* Neutralization of CXCL12 attenuates established pulmonary hypertension in rats. *Cardiovasc Res.* (2019).

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, 12/05/2022

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