

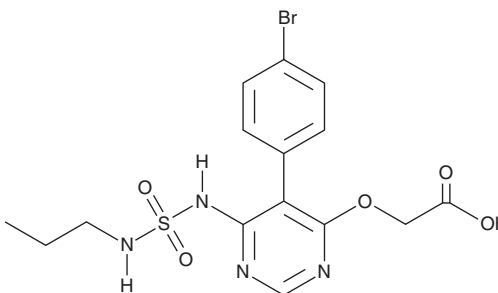
PRODUCT INFORMATION



ACT-373898

Item No. 34916

CAS Registry No.: 1433875-14-9
Formal Name: 2-[[5-(4-bromophenyl)-6-[[[(propylamino)sulfonyl]amino]-4-pyrimidinyl]oxy]-acetic acid
MF: C₁₅H₁₇BrN₄O₅S
FW: 445.3
Purity: ≥90%
UV/Vis.: λ_{max}: 257 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

ACT-373898 is supplied as a solid. A stock solution may be made by dissolving the ACT-373898 in the solvent of choice, which should be purged with an inert gas. ACT-373898 is soluble in ethanol and DMSO.

Description

ACT-373898 is a metabolite of the endothelin (ET) receptor type A (ET_A) and ET_B dual antagonist macitentan (Item No. 23304).¹

Reference

1. Dingemans, J., Sidharta, P.N., Maddrey, W.C., *et al.* Efficacy, safety and clinical pharmacology of macitentan in comparison to other endothelin receptor antagonists in the treatment of pulmonary arterial hypertension. *Expert Opin. Drug Saf.* **13**(3), 391-405 (2013).

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, 12/08/2022

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