

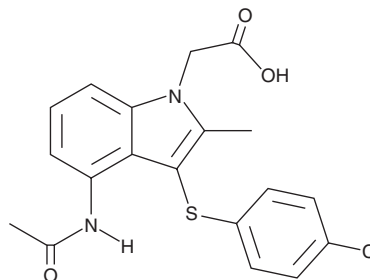
PRODUCT INFORMATION



AZD 1981

Item No. 20763

CAS Registry No.: 802904-66-1
Formal Name: 4-(acetylamino)-3-[(4-chlorophenyl)thio]-2-methyl-1H-indole-1-acetic acid
MF: C₁₉H₁₇ClN₂O₃S
FW: 388.9
UV/Vis.: λ_{max}: 223, 291 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

AZD 1981 is supplied as a crystalline solid. A stock solution may be made by dissolving the AZD 1981 in the solvent of choice, which should be purged with an inert gas. AZD 1981 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of AZD 1981 in these solvents is approximately 15 mg/ml.

AZD 1981 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AZD 1981 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. AZD 1981 has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

AZD 1981 is a DP₂/CRTH₂ receptor antagonist (IC₅₀ = 3.98 nM).¹ It is selective for DP₂ over a panel of receptors, ion channels, transporters, and enzymes, including DP₁, COX-1, and COX-2, at 10 μM. AZD 1981 (3-100 nM) inhibits increases in the expression of CD11b, a cell adhesion molecule, induced by 13,14-dihydro-15-keto-prostaglandin D₂ (13,14-dihydro-15-keto-PGD₂; Item No. 12610) in isolated human eosinophils. It also inhibits 13,14-dihydro-15-keto-PGD₂-induced chemotaxis of isolated human eosinophils and T cells in a concentration-dependent manner.

Reference

- Schmidt, J.A., Bell, F.M., Akam, E., *et al.* Biochemical and pharmacological characterization of AZD1981, an orally available selective DP₂ antagonist in clinical development for asthma. *Br. J. Pharmacol.* **168**(7), 1626-1638 (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, 12/19/2022

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