

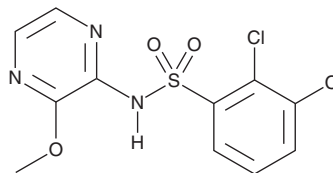
# PRODUCT INFORMATION



## AZD 2098

Item No. 27633

**CAS Registry No.:** 566203-88-1  
**Formal Name:** 2,3-dichloro-N-(3-methoxy-2-pyrazinyl)-benzenesulfonamide  
**MF:** C<sub>11</sub>H<sub>9</sub>Cl<sub>2</sub>N<sub>3</sub>O<sub>3</sub>S  
**FW:** 334.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 222 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 2098 is supplied as a crystalline solid. A stock solution may be made by dissolving the AZD 2098 in the solvent of choice. AZD 2098 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of AZD 2098 in these solvents is approximately 30 mg/ml.

AZD 2098 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AZD 2098 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. AZD 2098 has a solubility of approximately 0.20 mg/ml in a 1:4 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 2098 is a CC chemokine receptor 4 (CCR4) antagonist.<sup>1</sup> AZD 2098 inhibits CCL22-induced calcium flux in CHO cells expressing human recombinant CCR4 (IC<sub>50</sub> = 31.6 nM) and CCL22- and CCL17-mediated Th2 cell chemotaxis (IC<sub>50</sub> = 501.2 nM for both). AZD 2098 is selective for CCR4 over CXCR1, CXCR2, CCR1, CCR2b, CCR5, CCR7, and CCR8 up to 10 μM *in vitro*. Administration of AZD 2098 (0.07-2.5 mg/kg, p.o.) to ovalbumin-sensitized, antigen-challenged rats reduces alveolitis and leukocyte infiltration in the lung in a dose-dependent manner.

### Reference

1. Kindon, N.D., Andrews, G., Baxter, A., *et al.* Discovery of AZD-2098 and AZD-1678, two potent and bioavailable CCR4 receptor antagonists. *Med. Chem.* **8(9)**, 981-986 (2017).

#### 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, 10/04/2022

#### 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