

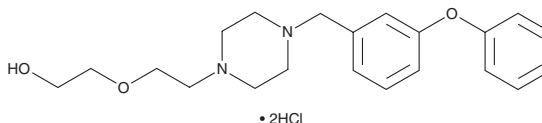
# PRODUCT INFORMATION



## ZK 756326 (hydrochloride)

Item No. 28454

**CAS Registry No.:** 1780259-94-0  
**Formal Name:** 2-[2-[4-[(3-phenoxyphenyl)methyl]-1-piperazinyl]ethoxy]-ethanol dihydrochloride  
**MF:**  $C_{21}H_{28}N_2O_3 \cdot 2HCl$   
**FW:** 429.4  
**Purity:**  $\geq 98\%$   
**Supplied as:** A crystalline solid  
**Storage:**  $-20^{\circ}C$   
**Stability:**  $\geq 4$  years



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

### Laboratory Procedures

ZK 756326 (hydrochloride) is supplied as a crystalline solid. Aqueous solutions of ZK 756326 (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of ZK 756326 (hydrochloride) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

ZK 756326 is a non-peptide agonist of the CC chemokine receptor CCR8. It inhibits binding of CCL1 to human CCR8 ( $IC_{50} = 1.8 \mu M$ ) and induces  $G_{ai}$ -mediated calcium-mobilization in U87 cells expressing the receptor.<sup>1</sup> It exhibits greater than 28-fold selectivity for CCR8 over 26 other GPCRs but also binds to the serotonin (5-HT) receptor subtypes 5-HT<sub>1A</sub>, 5-HT<sub>2B</sub>, 5-HT<sub>2C</sub>, and 5-HT<sub>5A</sub> with  $IC_{50}$  values of 5.4, 4.4, 34.8, and 16  $\mu M$ , respectively. ZK 756326 stimulates extracellular acidification in cells expressing human and mouse CCR8 ( $EC_{50}$ s = 254 nM and 2.6  $\mu M$ , respectively). ZK 756326 induces ERK phosphorylation, exhibits leukocyte chemotactic activity, and inhibits HIV fusion of cells expressing CD4 and CCR8.

### Reference

1. Haskell, C.A., Horuk, R., Liang, M., *et al.* Identification and characterization of a potent, selective nonpeptide agonist of the CC chemokine receptor CCR8. *Mol. Pharmacol.* **69**(1), 309-316 (2005).

#### 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, 11/10/2022

#### 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](http://WWW.CAYMANCHEM.COM)