

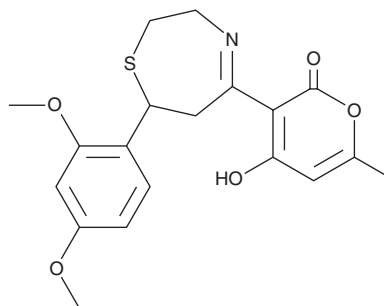
PRODUCT INFORMATION



KF38789

Item No. 43182

CAS Registry No.: 257292-29-8
Formal Name: 3-[7-(2,4-dimethoxyphenyl)-2,3,6,7-tetrahydro-1,4-thiazepin-5-yl]-4-hydroxy-6-methyl-2H-pyran-2-one
MF: C₁₉H₂₁NO₅S
FW: 375.4
Purity: ≥98%
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

KF38789 is supplied as a solid. A stock solution may be made by dissolving the KF38789 in the solvent of choice, which should be purged with an inert gas. KF38789 is soluble (≥10 mg/ml) in acetonitrile.

Description

KF38789 is an inhibitor of P-selectin-mediated cell adhesion.¹ It inhibits the binding of U937 cells to immobilized P-selectin immunoglobulin (IC₅₀ = 1.97 μM). It is selective for P-selectin-mediated cell adhesion over E- and or L-selectin-mediated adhesion of U937 and HL-60 cells, respectively, at 100 μM. KF38789 (0.1-10 μM) decreases P-selectin-induced production of superoxide by isolated human polymorphonuclear leukocytes (PMNs). It reduces the proliferation of T47D triple-negative breast and HCT116 colon cancer cells (EC₅₀s = 0.41 and 0.56 μM, respectively).² KF38789 (1 mg/kg) also reduces PMN infiltration into the peritoneum in a mouse model of thioglycollate-induced peritonitis.¹

References

1. Ohta, S., Inujima, Y., Abe, M., *et al.* Inhibition of P-selectin specific cell adhesion by a low molecular weight, non-carbohydrate compound, KF38789. *Inflamm. Res.* **50(11)**, 544-551 (2001).
2. Drewe, J., Kasibhatla, S., Tseng, B., *et al.* Discovery of 5-(4-hydroxy-6-methyl-2-oxo-2H-pyran-3-yl)-7-phenyl-(E)-2,3,6,7-tetrahydro-1,4-thiazepines as a new series of apoptosis inducers using a cell- and caspase-based HTS assay. *Bioorg. Med. Chem. Lett.* **17(17)**, 4987-4990 (2007).

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, 05/06/2025

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