

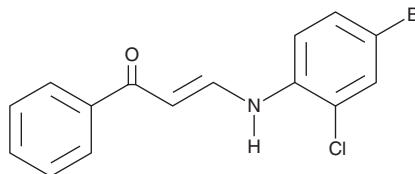
PRODUCT INFORMATION



SMI-6860766

Item No. 42960

CAS Registry No.: 433234-16-3
Formal Name: 3-[(4-bromo-2-chlorophenyl)amino]-1-phenyl-2-propen-1-one
MF: C₁₅H₁₁BrClNO
FW: 336.6
Purity: ≥98%
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

SMI-6860766 is supplied as a crystalline solid. A stock solution may be made by dissolving the SMI-6860766 in the solvent of choice, which should be purged with an inert gas. SMI-6860766 is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

SMI-6860766 is an inhibitor of the protein-protein interaction between TNF receptor-associated factor 6 (TRAF6) and CD40.¹ It binds to TRAF1, TRAF2, TRAF3, and TRAF6 (K_d s = 52, 30, 37, and 59 μ M, respectively). SMI-6860766 inhibits LPS-induced increases in the levels of NF- κ B in RAW 264.7 macrophages (IC_{50} = 0.3 μ M). It inhibits anti-CD40 antibody-induced increases in the levels of IL-1 β and IL-6 in primary mouse bone marrow-derived macrophages (BMDMs; IC_{50} s = 5.7 and 2.4 μ M, respectively). SMI-6860766 (10 μ mol/kg) increases survival in a mouse model of sepsis induced by cecal ligation and puncture. It decreases blood levels of glucose and adipose tissue levels of leukocytes, T helper cells (Ths), cytotoxic T cells, and M1 and M2 macrophages in a mouse model of diet-induced obesity (DIO) when administered at a dose of 10 μ mol/kg per day.²

References

1. Zarzycka, B., Seijkens, T., Nabuurs, S.B., et al. Discovery of small molecule CD40-TRAF6 inhibitors. *J. Chem. Inf. Model* **55**(2), 294-307 (2015).
2. van den Berg, S.M., Seijkens, T.T., Kusters, P.J., et al. Blocking CD40-TRAF6 interactions by small-molecule inhibitor 6860766 ameliorates the complications of diet-induced obesity in mice. *Int. J. Obes. (Lond)* **39**(5), 782-790 (2014).

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, 02/21/2025

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