

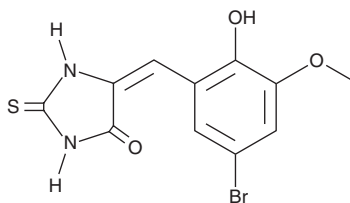
PRODUCT INFORMATION



(E/Z)-IT-603

Item No. 37792

CAS Registry No.: 292168-90-2
Formal Name: 5-[(5-bromo-2-hydroxy-3-methoxyphenyl)methylene]-2-thioxo-4-imidazolidinone
MF: $C_{11}H_9BrN_2O_3S$
FW: 329.2
Purity: $\geq 98\%$
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

(E/Z)-IT-603 is supplied as a solid. A stock solution may be made by dissolving the (E/Z)-IT-603 in the solvent of choice, which should be purged with an inert gas. (E/Z)-IT-603 is soluble in DMSO.

Description

(E/Z)-IT-603 is a mixture of (E)-IT-603 and (Z)-IT-603.¹ (Z)-IT-603 is a c-Rel inhibitor ($IC_{50} = 3 \mu\text{M}$). (Z)-IT-603 inhibits c-Rel DNA binding in an electrophoretic mobility-shift assay (EMSA). Preincubation of (Z)-IT-603 ($20 \mu\text{M}$) with mouse T cells prevents the development of graft-versus-host disease (GVHD) but does not affect graft-versus-tumor (GVT) activity in a mouse model of allogeneic hematopoietic stem cell transplantation.

References

1. Shono, Y., Tuckett, A.Z., Ouk, S., *et al.* A small-molecule c-Rel inhibitor reduces alloactivation of T cells without compromising antitumor activity. *Cancer Discov.* **4**(5), 578-591 (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, 06/17/2024

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