

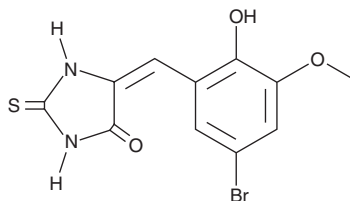
# PRODUCT INFORMATION



## (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<sub>11</sub>H<sub>9</sub>BrN<sub>2</sub>O<sub>3</sub>S  
**FW:** 329.2  
**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

(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 c-Rel inhibitor (IC<sub>50</sub> = 3 μM).<sup>1</sup> It inhibits c-Rel DNA binding to the *IL2* promoter in an EMSA. (E/Z)-IT-603 reduces the proliferation of OCI-Ly3 diffuse large B cell lymphoma cells (IC<sub>50</sub> = 18 μM). It reduces Huh7 epithelial carcinoma spheroid growth when used at a concentration of 20 μM in combination with the anthracycline antitumor antibiotic doxorubicin (Item No. 15007).<sup>2</sup> Preincubation of (E/Z)-IT-603 (20 μM) with murine T cells co-administered with A20 murine lymphoma cells to lethally irradiated mice as a model of T cell-depleted bone marrow transplantation reduces allo-stimulation without affecting anticancer T cell responses and increases survival.<sup>1</sup> *In vivo*, (E/Z)-IT-603 (24 mg/kg every other day), when used in combination with doxorubicin, reduces hepatic tumor burden, increases CD8<sup>+</sup> T cell and F4/80<sup>+</sup> macrophage tumor infiltrates, and increases tumoral mRNA encoding chemokine (C-X-C motif) ligand 9 (Cxcl9) in a Hep-53.4 murine hepatocellular carcinoma model.<sup>2</sup>

### References

- 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).
- Leslie, J., Hunter, J.E., Collins, A., *et al.* c-Rel-dependent Chk2 signaling regulates the DNA damage response limiting hepatocarcinogenesis. *Hepatology* **78**(4), 1050-1063 (2023).

#### 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, 10/30/2023

#### 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)