

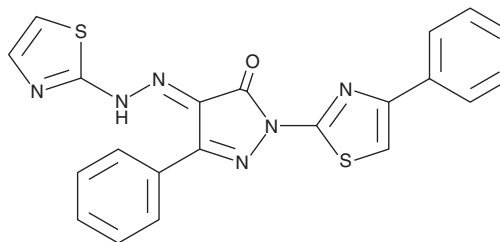
PRODUCT INFORMATION



BTSA1

Item No. 37305

CAS Registry No.: 314761-14-3
Formal Name: 3-phenyl-1-(4-phenyl-2-thiazolyl)-1H-pyrazole-4,5-dione, 4-[2-(2-thiazolyl)hydrazone]
Synonym: Bax Trigger Site Activator 1
MF: C₂₁H₁₄N₆OS₂
FW: 430.5
Purity: ≥98%
UV/Vis.: λ_{max}: 249 nm
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

BTSA1 is supplied as a solid. A stock solution may be made by dissolving the BTSA1 in the solvent of choice, which should be purged with an inert gas. BTSA1 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of BTSA1 in DMSO and DMF is approximately 2 and 10 mg/ml, respectively. BTSA1 is slightly soluble in ethanol.

Description

BTSA1 is an activator of Bax.¹ It binds to the Bax N-terminal activation site, also known as the trigger site (IC₅₀ = 250 nM), and induces Bax translocation membrane permeabilization in liposomes when used at concentrations ranging from 100 to 400 nM. BTSA1 (0.1-1 μM) induces apoptosis in various human and mouse acute myeloid leukemia (AML) cell lines. *In vivo*, BTSA (10 mg/kg) reduces bone marrow infiltration and promotes apoptosis in bone marrow infiltrates in a MOLM-13 AML mouse xenograft model.

Reference

1. Reyna, D.E., Garner, T.P., Lopez, A., *et al.* Direct activation of BAX by BTSA1 overcomes apoptosis resistance in acute myeloid leukemia. *Cancer Cell* **32**(4), 490-505 (2017).

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, 10/03/2023

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