

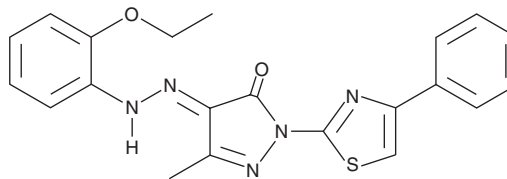
# PRODUCT INFORMATION



## BAM7

Item No. 14566

**CAS Registry No.:** 331244-89-4  
**Formal Name:** 3-methyl-1-(4-phenyl-2-thiazolyl)-1H-pyrazole-4,5-dione  
**Synonym:** Bax Activator Molecule 7  
**MF:** C<sub>21</sub>H<sub>19</sub>N<sub>5</sub>O<sub>2</sub>S  
**FW:** 405.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 236, 263, 426 nm  
**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

BAM7 is supplied as a crystalline solid. A stock solution may be made by dissolving the BAM7 in the solvent of choice, which should be purged with an inert gas. BAM7 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of BAM7 in these solvents is approximately 2 and 5 mg/ml, respectively.

BAM7 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BAM7 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. BAM7 has a solubility of approximately 0.33 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

BAM7 is a direct activator of Bax (EC<sub>50</sub> = 3.3 μM) that engages the Bax trigger site and promotes the functional oligomerization of Bax.<sup>1</sup> It is selective for the BH3 binding site on Bax and does not interact with BH3 binding pockets on other proteins. BAM7 induces Bax-dependent cell death in mouse embryo fibroblasts.<sup>1</sup>

### Reference

1. Gavathiotis, E., Reyna, D.E., Bellairs, J.A., *et al.* Direct and selective small-molecule activation of proapoptotic BAX. *Nat. Chem. Biol.* **8**(7), 639-645 (2012).

#### 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, 10/12/2022

#### 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