

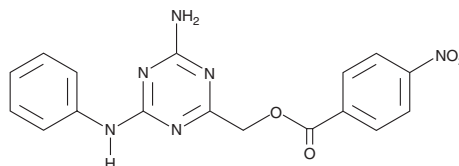
# PRODUCT INFORMATION



**TZ9**

Item No. 15964

**CAS Registry No.:** 1002789-86-7  
**Formal Name:** 4-amino-6-(phenylamino)-2-(4-nitrobenzoate)-1,3,5-triazine-2-methanol  
**MF:** C<sub>17</sub>H<sub>14</sub>N<sub>6</sub>O<sub>4</sub>  
**FW:** 366.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 209, 260 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

TZ9 is supplied as a crystalline solid. A stock solution may be made by dissolving the TZ9 in the solvent of choice. TZ9 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of TZ9 in these solvents is approximately 30 mg/ml.

TZ9 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, TZ9 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. TZ9 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

TZ9 is a cell-permeable triazine compound that inhibits the human E2 ubiquitin-conjugating enzyme Rad6B.<sup>1</sup> TZ9 selectively docks to the Rad6B catalytic site, inhibits Rad6B-induced histone H2A ubiquitination, downregulates intracellular β-catenin, induces G<sub>2</sub>/M arrest and apoptosis, and inhibits proliferation and migration of metastatic human breast cancer cells (IC<sub>50</sub> = 6 μM).<sup>1</sup>

## Reference

1. Sanders, M.A., Brahemi, G., Nangia-Makker, P., *et al.* Novel inhibitors of Rad6 ubiquitin conjugating enzyme: Design, synthesis, identification, and functional characterization. *Mol. Cancer Ther.* **12**(4), 373-383 (2013).

### 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, 11/21/2022

## 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](http://WWW.CAYMANCHEM.COM)