

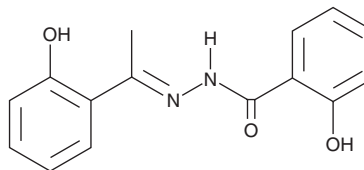
# PRODUCT INFORMATION



## Compound C108

Item No. 9002951

**CAS Registry No.:** 15533-09-2  
**Formal Name:** 2-hydroxy-benzoic acid, 2-[1-(2-hydroxyphenyl) ethylidene]hydrazide  
**MF:** C<sub>15</sub>H<sub>14</sub>N<sub>2</sub>O<sub>3</sub>  
**FW:** 270.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 286, 330 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Compound C108 is supplied as a crystalline solid. A stock solution may be made by dissolving the compound C108 in the solvent of choice. Compound C108 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of compound C108 in these solvents is approximately 1 mg/ml.

Compound C108 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, compound C108 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Compound C108 has a solubility of approximately 0.14 mg/ml in a 1:6 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Compound C108 is an inhibitor of the GTPase-activating protein (SH3 domain)-binding protein 2 (G3BP2), which is involved in breast tumor initiation.<sup>1</sup> In a mouse xenograft model of breast cancer, compound C108 (1 μM for 24 hours) reduces the proportion of tumor-initiating cells approximately 10-fold compared with untreated cells. In MDA-MB-453 cells, and in the patient-derived xenograft cell line MAXF 401, it diminishes mammosphere formation.

### Reference

1. Gupta, N., Badeaux, M., Liu, Y., *et al.* Stress granule-associated protein G3BP2 regulates breast tumor initiation. *Proc. Natl. Acad. Sci. USA* **114**(5), 1033-1038 (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, 06/14/2017

#### 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