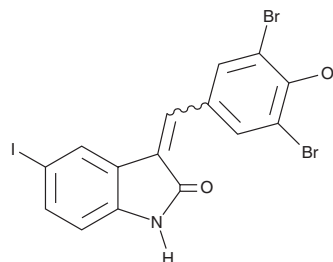


# PRODUCT INFORMATION

## GW 5074

Item No. 10010368

**CAS Registry No.:** 220904-83-6  
**Formal Name:** 3-[(3,5-dibromo-4-hydroxyphenyl)methylene]-1,3-dihydro-5-iodo-2H-indol-2-one  
**MF:** C<sub>15</sub>H<sub>8</sub>Br<sub>2</sub>INO<sub>2</sub>  
**FW:** 520.9  
**Purity:** ≥98% (mixture of *cis* and *trans*)  
**UV/Vis.:** λ<sub>max</sub>: 260, 348 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

GW 5074 is supplied as a crystalline solid. A stock solution may be made by dissolving the GW 5074 in the solvent of choice, which should be purged with an inert gas. GW 5074 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of GW 5074 in these solvents is approximately 15 mg/ml.

GW 5074 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GW 5074 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. GW 5074 has a solubility of approximately 0.1 mg/ml in a 1:8 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Raf-1 is a proto-oncogene serine/threonine protein kinase that signals from Ras to the MAPK/ERK signaling pathway.<sup>1-2</sup> This pathway mediates basic cellular functions, including proliferation, differentiation, and survival.<sup>1</sup> GW 5074 is a potent, selective, and cell-permeable inhibitor of Raf-1 (IC<sub>50</sub> = 9 nM).<sup>3</sup> It blocks phosphorylation of ERK1/2 by 90% in cells stimulated with epidermal growth factor when given at 5 μM.<sup>3</sup> GW 5074 shows more than 100-fold selectivity for Raf-1 versus several related kinases.<sup>3</sup>

### References

1. Dhillon, A.S. and Kolch, W. Untying the regulation of the Raf-1 kinase. *Arch. Biochem. Biophys.* **404**, 3-9 (2002).
2. Kunnimalaiyaan, M. and Chen, H. The Raf-1 pathway: a molecular target for treatment of select neuroendocrine tumors? *Anticancer Drugs* **17**, 139-142 (2006).
3. Lackey, K., Cory, M., David, R., et al. The discovery of potent cRaf1 kinase inhibitors. *Bioorg. Med. Chem. Lett.* **10**(3), 223-226 (2000).

#### 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, 12/05/2022

#### 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