

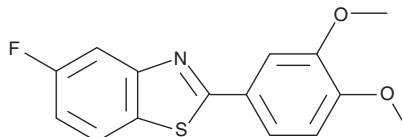
# Product Information



## GW 610

Catalog No. 10008313

**CAS Registry No.:** 872726-44-8  
**Formal Name:** 2-(3,4-dimethoxyphenyl)-5-fluorobenzothiazole  
**Synonym:** NSC 721648  
**MF:** C<sub>15</sub>H<sub>12</sub>FNO<sub>2</sub>S  
**FW:** 289.3  
**Purity:** ≥98%  
**Stability:** ≥2 years at -20°C  
**Supplied as:** A crystalline solid



### Laboratory Procedures

For long term storage, we suggest that GW 610 be stored as supplied at -20°C. It should be stable for at least two years.

GW 610 is supplied as a crystalline solid. A stock solution may be made by dissolving the GW 610 in an organic solvent purged with an inert gas. GW 610 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of GW 610 in DMSO is approximately 0.5 mg/ml and 10 mg/ml in DMF.

GW 610 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GW 610 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. GW 610 has a solubility of approximately 0.25 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.

GW 610 is an antitumor benzothiazole that shows growth-inhibitory activity against several cancer cell lines. In MCF-7 and MDA 468 human cancer cell lines, potent antiproliferative activity (growth inhibition (GI<sub>50</sub>) < 0.1 nM) was observed.<sup>1</sup> Potent and selective activity was also observed in the NCI 60 human cancer cell line panel. In contrast with previously reported 2-(4-aminophenyl)benzothiazoles, GW 610 is not reliant on induction of cytochrome P 1A1 (CYP1A1) expression for antitumor activity.

### Reference

1. Mortimer, C.G., Wells, G., Crochard, J.-P., *et al.* Antitumor benzothiazoles. 26. 2-(3,4-Dimethoxyphenyl)-5-fluorobenzothiazole (GW 610, NSC 721648), a simple fluorinated 2-arylbenzothiazole, shows potent and selective inhibitory activity against lung, colon, and breast cancer cell lines. *J. Med. Chem.* **49**, 179-185 (2006).

### Related Products

PRIMA-1 - Cat. No. 63520 • Nutlin-3 - Cat. No. 10004372 • RITA - Cat. No. 10006426 • OSU03012 - Cat. No. 10008005

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY; NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 05/07/2010

### Cayman Chemical

#### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

#### Phone

(800) 364-9897  
(734) 971-3335

#### Fax

(734) 971-3640

#### E-Mail

custserv@caymanchem.com

#### Web

www.caymanchem.com