

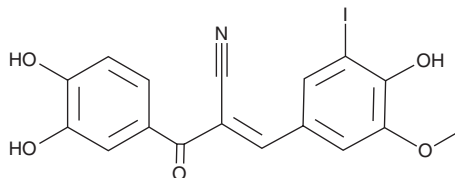
# PRODUCT INFORMATION



## I-OMe AG-538

Item No. 35090

**CAS Registry No.:** 1094048-77-7  
**Formal Name:** 3,4-dihydroxy- $\alpha$ E-[(4-hydroxy-3-iodo-5-methoxyphenyl)methylene]- $\beta$ -oxo-benzenepropanenitrile  
**Synonym:** I-OMe Tyrphostin AG-538  
**MF:** C<sub>17</sub>H<sub>12</sub>INO<sub>5</sub>  
**FW:** 437.2  
**Purity:**  $\geq 95\%$   
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:**  $\geq 4$  years



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

### Laboratory Procedures

I-OMe AG-538 is supplied as a solid. A stock solution may be made by dissolving the I-OMe AG-538 in the solvent of choice, which should be purged with an inert gas. I-OMe AG-538 is soluble ( $\geq 10$  mg/ml) in DMSO.

### Description

I-OMe AG-538 is an inhibitor of the insulin-like growth factor 1 receptor (IGF-1R;  $IC_{50} = 3.4 \mu M$ ) and phosphatidylinositol 5-phosphate 4-kinase  $\alpha$  (PI5P4K $\alpha$ ;  $IC_{50} = 1 \mu M$ ).<sup>1,2</sup> It is selectively cytotoxic to nutrient-deprived PANC-1 pancreatic cancer cells over nutrient-sufficient PANC-1 cells.<sup>3</sup>

### References

1. Blum, G., Gazit, A., and Levitzki, A. Substrate competitive inhibitors of IGF-1 receptor kinase. *Biochemistry* **39**(51), 15705-15712 (2000).
2. Davis, M.I., Sasaki, A.T., Shen, M., *et al.* A homogeneous, high-throughput assay for phosphatidylinositol 5-phosphate 4-kinase with a novel, rapid substrate preparation. *PLoS One* **8**(1), e54127 (2013).
3. Momose, I., Kunitomo, S., Osono, M., *et al.* Inhibitors of insulin-like growth factor-1 receptor tyrosine kinase are preferentially cytotoxic to nutrient-deprived pancreatic cancer cells. *Biochem. Biophys. Res. Commun.* **380**(1), 171-176 (2009).

#### 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, 12/05/2024

#### 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)