

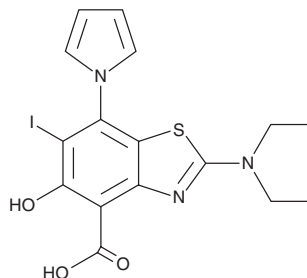
PRODUCT INFORMATION



MB710

Item No. 45499

CAS Registry No.: 2230044-57-0
Formal Name: 2-(diethylamino)-5-hydroxy-6-iodo-7-(1H-pyrrol-1-yl)-4-benzothiazolecarboxylic acid
MF: C₁₆H₁₆IN₃O₃S
FW: 457.3
Purity: ≥98%
Supplied as: A 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

MB710 is supplied as a solid. A stock solution may be made by dissolving the MB710 in the solvent of choice, which should be purged with an inert gas. MB710 is sparingly soluble (1-10 mg/ml) in acetonitrile.

Description

MB710 is a stabilizer of mutant p53.¹ It increases the thermal stability and inhibits aggregation of the mutant p53^{Y220C} DNA-binding domain ($K_d = 4 \mu\text{M}$). MB710 selectively inhibits growth of NUGC-3 gastric adenocarcinoma cells expressing p53^{Y220C} over WI38 cells that do not express the p53^{Y220C} mutant ($IC_{50s} = 90$ and $>120 \mu\text{M}$, respectively). It also selectively reduces the viability of Huh7 p53^{Y220C}-expressing human epithelial carcinoma cells over Huh7 p53^{-/-} cells at 90 and 120 μM , respectively.

Reference

1. Baud, M.G.J., Bauer, M.R., Verduci, L., *et al.* Aminobenzothiazole derivatives stabilize the thermolabile p53 cancer mutant Y220C and show anticancer activity in p53-Y220C cell lines. *Eur. J. Med. Chem.* **152**, 101-114 (2018).

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, 04/03/2026

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