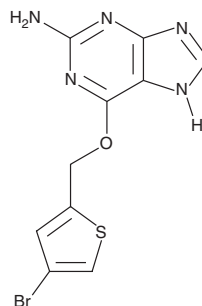


# PRODUCT INFORMATION



## Lomeguatrib Item No. 11732

**CAS Registry No.:** 192441-08-0  
**Formal Name:** 6-[(4-bromo-2-thienyl)methoxy]-  
9H-purin-2-amine  
**Synonym:** PaTrin 2  
**MF:** C<sub>10</sub>H<sub>8</sub>BrN<sub>5</sub>OS  
**FW:** 326.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 239, 284 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

Lomeguatrib is supplied as a crystalline solid. A stock solution may be made by dissolving the lomeguatrib in the solvent of choice. Lomeguatrib is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of lomeguatrib in these solvents is approximately 5 and 10 mg/ml, respectively.

Lomeguatrib is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

### Description

O<sup>6</sup>-Methylguanine-DNA methyltransferase (MGMT) (Item No. 11176) is a DNA repair protein which removes an alkyl group from the O<sup>6</sup> position on guanine in an autoinactivating reaction.<sup>1</sup> Although important in normal DNA repair, this reaction confers resistance to treatments that use O<sup>6</sup>-alkylating agents to produce cytotoxicity, e.g., in cancer.<sup>2,3</sup> Lomeguatrib is a modified guanine base which acts as a pseudosubstrate inactivator of MGMT (IC<sub>50</sub> = ~3 nM).<sup>1,2</sup> A non-toxic compound, lomeguatrib completely inactivates MGMT in human prostate and colorectal tumors when given as a single 120 mg oral dose and in primary central nervous system cancers at 160 mg.<sup>4,5</sup>

### References

1. McMurry, T.B.H. MGMT inhibitors-the Trinity College-Paterson Institute experience, a chemist's perception. *DNA Repair* **6(8)**, 1161-1169 (2007).
2. Clemons, M., Kelly, J., Watson, A.J., et al. O<sup>6</sup>-(4-bromothienyl)guanine reverses temozolomide resistance in human breast tumour MCF-7 cells and xenografts. *Br. J. Cancer* **93**, 1152-1156 (2005).
3. Marchesi, F., Turriziani, M., Tortorelli, G., et al. Triazene compounds: Mechanism of action and related DNA repair systems. *Pharmacol. Res.* **56(4)**, 275-287 (2007).
4. Ranson, M., Middleton, M.R., Bridgewater, J., et al. Lomeguatrib, a potent inhibitor of O<sup>6</sup>-alkylguanine-DNA-alkyltransferase: Phase I safety, pharmacodynamic, and pharmacokinetic trial and evaluation in combination with temozolomide in patients with advanced solid tumors. *Clin. Cancer Res.* **12**, 1577-1584 (2006).
5. Watson, A.J., Sabharwal, A., Thorncroft, M., et al. Tumor O<sup>6</sup>-methylguanine-DNA methyltransferase inactivation by oral lomeguatrib. *Clin. Cancer Res.* **16(2)**, 743-749 (2010).

#### 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, 11/08/2022

#### 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