

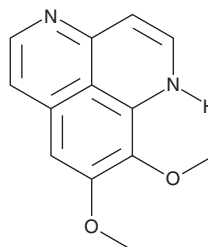
# PRODUCT INFORMATION



## Aptamine

Item No. 24297

**CAS Registry No.:** 85547-22-4  
**Formal Name:** 8,9-dimethoxy-1H-benzo[de][1,6]naphthyridine  
**MF:** C<sub>13</sub>H<sub>12</sub>N<sub>2</sub>O<sub>2</sub>  
**FW:** 228.3  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -80°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

Aptamine is supplied as a solid. A stock solution may be made by dissolving the aptamine in the solvent of choice. Aptamine is soluble in organic solvents such as ethanol and DMSO, which should be purged with an inert gas.

### Description

Aptamine is a marine sponge alkaloid originally isolated from *A. aptos* with diverse biological activities, including antiproliferative and antidepressant properties.<sup>1-4</sup> It is an agonist of  $\delta$ - and  $\mu$ -opioid receptors (EC<sub>50</sub>s = 5.1 and 10.1  $\mu$ M, respectively, in HEK293 cells expressing human recombinant receptors) and a competitive antagonist of  $\alpha$ -adrenergic receptors ( $\alpha$ -ARs; pA<sub>2</sub>s = 4.88 and 5.43, respectively, in isolated rabbit aorta and renal artery).<sup>1,4</sup> Aptamine inhibits growth of HeLa cervical cancer (IC<sub>50</sub> = 15  $\mu$ g/ml) and K562 leukemia (GI<sub>50</sub> = 10  $\mu$ M) cells and induces cell cycle arrest in the G<sub>2</sub>/M phase in K562 leukemia and MG63 osteosarcoma cells.<sup>2,3,5</sup> It is a proteasome inhibitor that inhibits chymotrypsin-, caspase-, and trypsin-like activity in a partially purified rat liver proteasome preparation (IC<sub>50</sub>s = 1.6, 2.7, and 18  $\mu$ g/ml, respectively) and dose-dependently activates p21 independent of p53 in MG63 cells when used at concentrations ranging from 20 to 50  $\mu$ M.<sup>2,5</sup> Aptamine (40 mg/kg, i.p.) decreases immobility time in the forced swim test in mice, indicating antidepressant-like activity.<sup>4</sup>

### References

- Ohizumi, Y., Kajiwarra, A., Nakamura, H., *et al.*  $\alpha$ -Adrenoceptor blocking action of aptamine, a novel marine natural product, in vascular smooth muscle. *J. Pharm. Pharmacol.* **36(11)**, 785-786 (1984).
- Tsukamoto, S., Yamanokuchi, R., Yoshitomi, M., *et al.* Aptamine, an alkaloid from the sponge *Aptos suberitoides*, functions as a proteasome inhibitor. *Bioorg. Med. Chem. Lett.* **20(11)**, 3341-3343 (2010).
- Jin, M., Zhao, W., Zhang, Y., *et al.* Antiproliferative effect of aptamine on human chronic myeloid leukemia K562 cells. *Int. J. Mol. Sci.* **12(11)**, 7352-7359 (2011).
- Johnson, T.A., Milan-Lobo, L., Che, T., *et al.* Identification of the first marine-derived opioid receptor "balanced" agonist with a signaling profile that resembles the endorphins. *ACS Chem. Neurosci.* **8(3)**, 473-485 (2017).
- Aoki, S., Kong, D., Suna, H., *et al.* Aptamine, a spongean alkaloid, activates p21 promoter in a p53-independent manner. *Biochem. Biophys. Res. Commun.* **342(1)**, 101-106 (2006).

#### 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, 12/13/2018

#### 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