

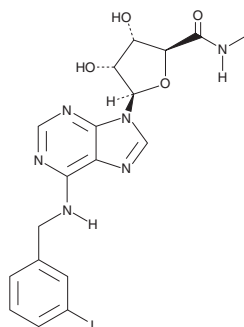
# PRODUCT INFORMATION



## Piclidenoson

Item No. 28445

**CAS Registry No.:** 152918-18-8  
**Formal Name:** 1-deoxy-1-[6-[[[3-iodophenyl]methyl]amino]-9H-purin-9-yl]-N-methyl-β-D-ribofuranuronamide  
**Synonyms:** CF-101, IB-MECA  
**MF:** C<sub>18</sub>H<sub>19</sub>IN<sub>6</sub>O<sub>4</sub>  
**FW:** 510.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 268 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

Piclidenoson is supplied as a crystalline solid. A stock solution may be made by dissolving the piclidenoson in the solvent of choice, which should be purged with an inert gas. Piclidenoson is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of piclidenoson in these solvents is approximately 2 mg/ml.

Piclidenoson is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, piclidenoson should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Piclidenoson has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Piclidenoson is an adenosine A<sub>3</sub> receptor agonist (K<sub>i</sub> = 1.7 nM).<sup>1</sup> It is selective for adenosine A<sub>3</sub> over A<sub>1</sub> and A<sub>2</sub> receptors (K<sub>i</sub>s = 55.3 and 3,099 nM, respectively). *In vivo*, piclidenoson (100 μg/kg) reduces wall thickening in a rabbit model of ischemia-reperfusion injury induced by coronary artery occlusion, an effect that can be reversed by the adenosine receptor antagonist 8-*p*-sulfophenyl theophylline.<sup>2</sup> It reduces hepatocellular tumor growth, liver inflammation, and neuropathy in a rat model of bone-residing breast cancer and decreases the number of lung metastases in a model of metastatic murine melanoma.<sup>1</sup> Piclidenoson (1 and 3 mg/kg) also reduces colonic inflammatory cell infiltration and damage in a mouse model of colitis induced by dextran sulfate (DSS; Item No. 23250) and in *IL-10*<sup>-/-</sup> mice.<sup>3</sup>

### References

1. Jacobson, K.A., Merighi, S., Varani, K., *et al.* A<sub>3</sub> adenosine receptors as modulators of inflammation: From medicinal chemistry to therapy. *Med. Res. Rev.* **38(4)**, 1031-1072 (2018).
2. Auchampach, J.A., Rizvi, A., Qiu, Y., *et al.* Selective activation of A<sub>3</sub> adenosine receptors with N<sup>6</sup>-(3-iodobenzyl)adenosine-5'-N-methyluronamide protects against myocardial stunning and infarction without hemodynamic changes in conscious rabbits. *Circ. Res.* **80(6)**, 800-809 (1997).
3. Mabley, J., Soriano, F., Pacher, P., *et al.* The adenosine A<sub>3</sub> receptor agonist, N<sup>6</sup>-(3-iodobenzyl)-adenosine-5'-N-methyluronamide, is protective in two murine models of colitis. *Eur. J. Pharmacol.* **466(3)**, 323-329 (2003).

#### 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, 06/08/2020

#### 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