

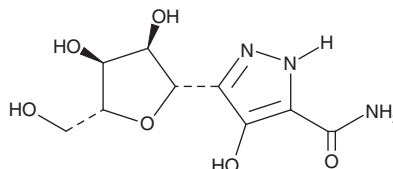
# PRODUCT INFORMATION



## Pyrazofurin

Item No. 43569

**CAS Registry No.:** 30868-30-5  
**Formal Name:** 4-hydroxy-3-β-D-ribofuranosyl-1H-pyrazole-5-carboxamide  
**MF:** C<sub>9</sub>H<sub>13</sub>N<sub>3</sub>O<sub>6</sub>  
**FW:** 259.2  
**Purity:** ≥95%  
**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

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

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of pyrazofurin can be prepared by directly dissolving the solid in aqueous buffers. Pyrazofurin is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

### Description

Pyrazofurin is a nucleoside derivative that has been found in *S. candidus* and has antiviral and anticancer activities.<sup>1-3</sup> It is an inhibitor of orotidine-5'-monophosphate decarboxylase (OMP decarboxylase) that inhibits DNA synthesis.<sup>2</sup> Pyrazofurin inhibits the cytopathic effect of parainfluenza-3 virus, measles virus, vaccinia virus, and herpes simplex virus 2 (HSV-2) in Vero cells (EC<sub>50</sub>s = 3, 9.6, 2.8, and 20 μM, respectively).<sup>1</sup> It also reduces infection of Calu-3 cells by severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) with a synergistic effect when used in combination with molnupiravir (Item No. 29586) or remdesivir (Item No. 30354).<sup>3</sup> Pyrazofurin reduces tumor growth in L1210 and P388 leukemia, Lewis lung, and B16 murine melanoma models (ID<sub>50</sub>s = 3.2, 3.3, 1.8, and 63 nM, respectively).<sup>1</sup>

### References

1. Petrie, C.R., 3rd., Revankar, G.R., Dalley, N.K., *et al.* Synthesis and biological activity of certain nucleoside and nucleotide derivatives of pyrazofurin. *J. Med. Chem.* **29**(2), 268-278 (1986).
2. Ringer, D.P., Howell, B.A., and Etheredge, J.L. Alteration in de novo pyrimidine biosynthesis during uridine reversal of pyrazofurin-inhibited DNA synthesis. *J. Biochem. Toxicol.* **69**(1), 19-27 (1991).
3. Schultz, D.C., Johnson, R.M., Ayyanathan, K., *et al.* Pyrimidine inhibitors synergize with nucleoside analogues to block SARS-CoV-2. *Nature* **604**(7904), 134-140 (2022).

#### 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, 09/08/2025

#### 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