

# 5-Iodotubercidin

Protein kinase inhibitor

Inhibits ERK2 (Ki = 525nM). Also inhibits adenosine kinase (Ki = 30nM), CK1 and CK2 and insulin receptor kinase (IC50 ranging from 0.4-28µM).

Citations: 2

[View Online »](#)

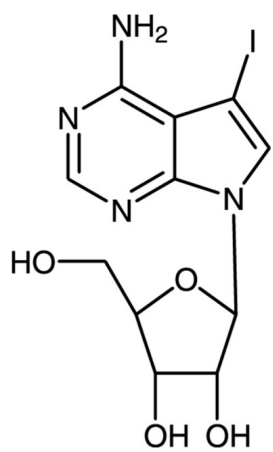
Ordering Information

[Order Online »](#)

BML-EI293-0001	1mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	4-Amino-5-iodo-7-(β-D-ribofuranosyl)-pyrrolo[2,3-d]pyrimidine
<b>Appearance</b>	Off-white solid
<b>CAS</b>	24386-93-4
<b>Couple Target</b>	Adenosine kinase, Casein kinase, Erk, Insulin receptor
<b>Couple Type</b>	Inhibitor
<b>Formula</b>	$C_{11}H_{13}INO_4$
<b>Identity</b>	Determined by 1H-NMR and MS
<b>MW</b>	392.1
<b>Purity</b>	≥98% (TLC)
<b>Solubility</b>	Soluble in DMSO (25mg/ml) or 100% ethanol (8mg/ml warm)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)