

# PRODUCT INFORMATION

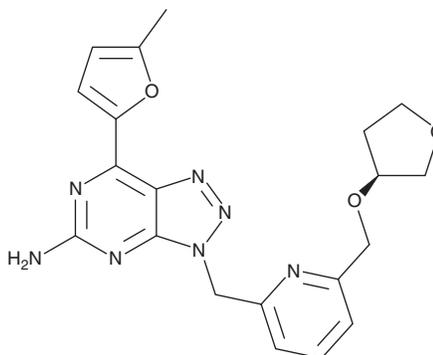


## CPI-444

Item No. 25024

**CAS Registry No.:** 1202402-40-1  
**Formal Name:** 7-(5-methyl-2-furanyl)-3-[[[6-[[[(3S)-tetrahydro-3-furanyl]oxy]methyl]-2-pyridinyl]methyl]-3H-1,2,3-triazolo[4,5-d]pyrimidin-5-amine

**Synonym:** Ciforadenant  
**MF:** C<sub>20</sub>H<sub>21</sub>N<sub>7</sub>O<sub>3</sub>  
**FW:** 407.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 321, 352 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

CPI-444 is supplied as a crystalline solid. A stock solution may be made by dissolving the CPI-444 in the solvent of choice, which should be purged with an inert gas. CPI-444 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CPI-444 in these solvents is approximately 30 mg/ml.

CPI-444 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CPI-444 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CPI-444 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

CPI-444 is an antagonist of the adenosine A<sub>2A</sub> receptor.<sup>1</sup> It reduces tumor area in mouse model of murine Her2/neu-expressing breast cancer.

### Reference

1. Scott, B.A., Armstrong, T., and Jaffee, E.M. Inhibition of adenosine A<sub>2A</sub> receptor (A<sub>2A</sub>R) by CPI-444 enhances CD8<sup>+</sup> T cell killing of a HER-2/neu expressing murine tumor. *Cancer Res.* **74(14)**, 320 (2016).

#### 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, 11/14/2022

#### 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