

# PRODUCT INFORMATION



## AM2394

Item No. 20992

**CAS Registry No.:** 1442684-77-6  
**Formal Name:** N-[6'-(2-hydroxy-2-methylpropoxy)-4-[(5-methyl-3-pyridinyl)oxy][3,3'-bipyridin]-6-yl]-N'-methyl-urea

**MF:** C<sub>22</sub>H<sub>25</sub>N<sub>5</sub>O<sub>4</sub>

**FW:** 423.5

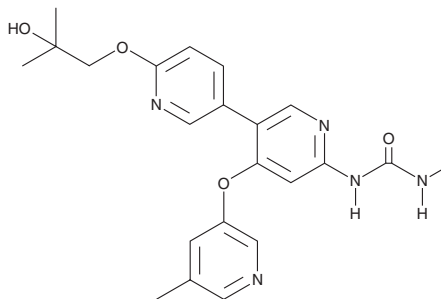
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 269 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

AM2394 is supplied as a crystalline solid. A stock solution may be made by dissolving the AM2394 in the solvent of choice, which should be purged with an inert gas. AM2394 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of AM2394 in these solvents is approximately 1 mg/ml.

AM2394 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AM2394 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. AM2394 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

AM2394 is an activator of glucokinase with an EC<sub>50</sub> value of 60 nM in an enzyme assay in the presence of glucose.<sup>1</sup> Oral administration of AM2394 (3-30 mg/kg) decreases plasma glucose levels in an oral glucose tolerance test in an *ob/ob* mouse model of diabetes.

### References

1. Dransfield, P.J., Pattaropong, V., Lai, S., *et al.* Novel series of potent glucokinase activators leading to the discovery of AM-2394. *ACS Med. Chem. Lett.* **7**(7), 714-718 (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, 10/10/2022

#### 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