

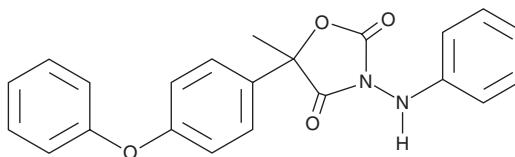
# PRODUCT INFORMATION



## Famoxadone

Item No. 34470

**CAS Registry No.:** 131807-57-3  
**Formal Name:** 5-methyl-5-(4-phenoxyphenyl)-3-(phenylamino)-2,4-oxazolidinedione  
**Synonym:** DPX-JE874  
**MF:** C<sub>22</sub>H<sub>18</sub>N<sub>2</sub>O<sub>4</sub>  
**FW:** 374.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 229 nm  
**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

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

Famoxadone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, famoxadone should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Famoxadone has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Famoxadone is a fungicide.<sup>1</sup> It inhibits mitochondrial complex III, also known as the cytochrome bc<sub>1</sub> complex, and electron transport in rat heart submitochondria (IC<sub>50</sub>s = 28 and 4.5 ng/ml, respectively). It reduces oxygen consumption in, and induces disintegration of, *P. infestans* zoospores when used at a concentration of 2 µg/ml. Famoxadone inhibits the growth of *P. viticola* on grape seedlings (EC<sub>50</sub> = 3 µg/ml). Formulations containing famoxadone have been used to control various fungal diseases in crops.

### Reference

1. Jordan, D.B., Livingston, R.S., Bisaha, J.J., *et al.* Mode of action of famoxadone. *Pestic. Sci.* 55, 105-118 (1999).

#### 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, 11/23/2022

#### 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