

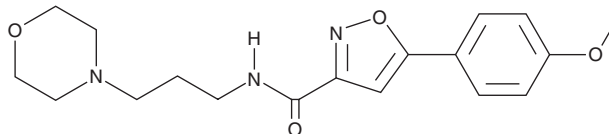
PRODUCT INFORMATION



CFTR Corrector 11

Item No. 45309

CAS Registry No.: 688050-45-5
Formal Name: 5-(4-methoxyphenyl)-N-[3-(4-morpholinyl)propyl]-3-isoxazolecarboxamide
MF: C₁₈H₂₃N₃O₄
FW: 345.4
Purity: ≥98%
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

CFTR corrector 11 is supplied as a solid. A stock solution may be made by dissolving the CFTR corrector 11 in the solvent of choice, which should be purged with an inert gas. CFTR corrector 11 is soluble (≥10 mg/ml) in DMSO. CFTR corrector 11 is sparingly soluble (1-10 mg/ml) in ethanol.

Description

CFTR corrector 11 is a substituted isoxazole containing morpholino and methoxy-benzene moieties.

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, 05/05/2026

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