

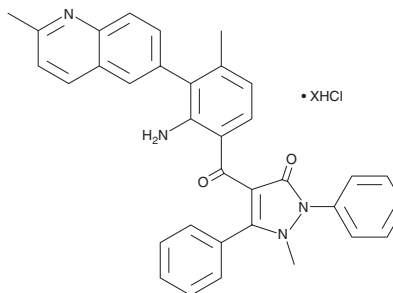
PRODUCT INFORMATION



AMG PERK 44 (hydrochloride)

Item No. 19126

Formal Name: 4-(2-amino-4-methyl-3-(2-methylquinolin-6-yl)benzoyl)-1-methyl-2,5-diphenyl-1,2-dihydro-3H-pyrazol-3-one, hydrochloride
MF: $C_{34}H_{28}N_4O_2 \cdot XHCl$
FW: 524.6
Purity: $\geq 98\%$
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

AMG PERK 44 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the AMG PERK 44 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. AMG PERK 44 (hydrochloride) is sparingly soluble (1-10 mg/ml) in methanol.

Description

AMG PERK 44 is a protein kinase R-like ER kinase (PERK) inhibitor ($IC_{50} = 6$ nM).¹ It is selective for PERK over general control nonderepressible 2 kinase (GCN2) and B-RAF ($IC_{50}s = 7,300$ and $>1,000$ nM, respectively). It inhibits PERK autophosphorylation in a cell-free assay ($IC_{50} = 84$ nM) and rescues protein synthesis in U2OS cells ($EC_{50} = 400$ nM). AMG PERK 44 (1, 3, 10, and 30 mg/kg) reduces intratumoral PERK autophosphorylation in a doxycycline-inducible PERK Matrigel™ plug assay.

Reference

1. Smith, A.L., Andrews, K.L., Beckmann, H., *et al.* Discovery of 1H-pyrazol-3(2H)-ones as potent and selective inhibitors of protein kinase R-like endoplasmic reticulum kinase (PERK). *J. Med. Chem.* **58**(3), 1426-1441 (2015).

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, 08/01/2025

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