

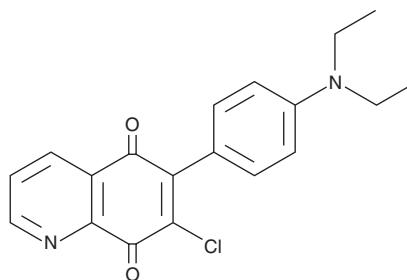
PRODUCT INFORMATION



EA4

Item No. 10111

CAS Registry No.: 389614-94-2
Formal Name: 7-chloro-6-[4-(diethylamino)phenyl]-5,8-quinolinedione
MF: C₁₉H₁₇ClN₂O₂
FW: 340.8
Purity: ≥95%
UV/Vis.: λ_{max}: 202, 242, 280, 600 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

EA4 is supplied as a crystalline solid. A stock solution may be made by dissolving the EA4 in the solvent of choice, which should be purged with an inert gas. EA4 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of EA4 in ethanol is approximately 1 mg/ml and approximately 20 mg/ml in DMSO and DMF.

EA4 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, EA4 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. EA4 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

rPLA₂ is a calcium-dependent cytosolic phospholipase A₂ (cPLA₂) initially isolated and characterized from bovine and human red blood cells (RBCs).¹ It has a molecular mass of 42 kDa and biochemical properties similar to cPLA₂ Type IV. EA4 is a quinone derivative that inhibits rPLA₂ with a K_i of 130 μM. It inhibits the ionophore-induced arachidonic acid release from human and bovine red blood cells (RBCs), indicating that rPLA₂ is responsible for the Ca²⁺-dependent release of arachidonic acid from mammalian RBCs.¹ EA4 causes significant time- and concentration-dependent induction of cytochrome P450 1A1 (CYP1A1) mRNA and protein in murine hepatoma Hepa-1c1c7 cells.² EA4 also modulates CYP1A1 and CYP1B1 expression in other cell lines such as MCF-7, Hep-G₂, and HL-60.²

References

1. Shin, H.S., Chin, M.R., Kim, J.S., *et al.* Purification and characterization of a cytosolic, 42-kDa and Ca²⁺-dependent phospholipase A₂ from bovine red blood cells. *J. Biol. Chem.* **277**(23), 21086-21094 (2002).
2. Chun, Y.J., Lee, B.Y., Yang, S.A., *et al.* Induction of cytochrome P450 1A1 gene expression by a vitamin K₃ analog in mouse hepatoma Hepa-1c1c7 cells. *Mol. Cells* **12**(2), 190-196 (2001).

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, 03/27/2023

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