

# PRODUCT INFORMATION



**CAY10648**

Item No. 9000549

**CAS Registry No.:** 916048-02-7  
**Formal Name:** 5-chloro-2,2'-dioxo-1,1-dimethylethyl ester-spiro[3H-indole-3,3'-pyrrolidine]-1(2H)-acetic acid

**MF:** C<sub>17</sub>H<sub>19</sub>ClN<sub>2</sub>O<sub>4</sub>

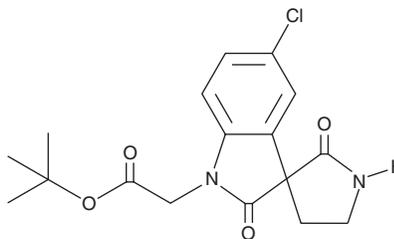
**FW:** 350.8

**Purity:** ≥98%

**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

CAY10648 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10648 in the solvent of choice, which should be purged with an inert gas. CAY10648 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of CAY10648 in these solvents is approximately 1, 10, and 5 mg/ml, respectively.

CAY10648 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10648 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CAY10648 has a solubility of approximately 0.25 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

CAY10648 is an intermediate used in the synthesis of potent DP<sub>2</sub> receptor antagonists based on a novel spiro-indolinone compound.<sup>1</sup>

## Reference

1. Crosignani, S., Page, P., Missotten, M., *et al.* Discovery of a new class of potent, selective, and orally bioavailable CRTH2 (DP2) receptor antagonists for the treatment of allergic inflammatory diseases. *J. Med. Chem.* **51**(7), 2227-2243 (2008).

### 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, 04/10/2024

## 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