

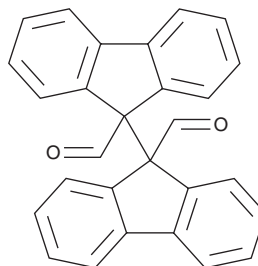
# PRODUCT INFORMATION



## Raptinal

Item No. 39131

**CAS Registry No.:** 1176-09-6  
**Formal Name:** [9,9'-Bi-9H-fluorene]-9,9'-dicarboxaldehyde  
**Synonyms:** 9H,9'H-9,9'-Bifluorene-9,9'-dicarbaldehyde, NSC 610970, SML1745  
**MF:** C<sub>28</sub>H<sub>18</sub>O<sub>2</sub>  
**FW:** 386.4  
**Purity:** ≥95%  
**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

Raptinal is supplied as a solid. A stock solution may be made by dissolving the raptinal in the solvent of choice, which should be purged with an inert gas. Raptinal is slightly soluble (0.1-1 mg/ml) in ethanol and DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of raptinal can be prepared by directly dissolving the solid in aqueous buffers. Raptinal is slightly soluble (0.1-1 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

### Description

Raptinal is an inducer of apoptosis.<sup>1</sup> It is hydrolyzed into its active form in aqueous solutions. Raptinal (10 µM) induces apoptosis in 17 cancerous and HFF-1 IRR, MCF-10A, MEFs, and 3T3-SA non-cancerous cells. *In vivo*, raptinal (20 mg/kg per day) decreases tumor mass in a 4T1 murine breast cancer model.

### Reference

1. Palchaudhuri, R., Lambrecht, M.J., Botham, R.C., *et al.* A small molecule that induces intrinsic pathway apoptosis with unparalleled speed. *Cell Rep.* **13**(9), 2027-2036 (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, 05/21/2024

#### 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