

# Prostaglandin A2

## CDK inhibitor

Blocks cell cycle at the G1 and G2/M phase [1] by inhibiting G1 phase cyclin dependent kinases [2] downregulating cyclin D1 [3]. Modulates all three aspects of the glutathione-mediated biotransformation system, i.e. GSH levels, glutathione S-transferase activity and transport of GSH conjugates [4]. Antitumor activity. Induces  $\gamma$ -glutamylcysteine synthetase and apoptosis in L-1210 cells.

Citations: 8

[View Online »](#)

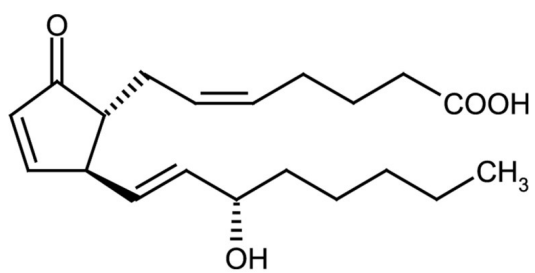
## Ordering Information

[Order Online »](#)

BML-PG002-0010	10mg
----------------	------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C.
Handling	Store at -20°C or colder. For experiments with cells or tissues, dissolve in ethanol and add aliquots directly to aqueous media. For HPLC, aliquots in acetone solution should be evaporated to dryness under a stream of nitrogen and HPLC mobile phase added to the desired concentration. Avoid basic conditions.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Medullin, Prosta-5,10,13-trien-1-oic acid, 15-hydroxy-9-oxo-, (5Z,13E,15S)-, PGA2
Appearance	Neat oil.
CAS	13345-50-1
Couple Target	CDK
Couple Type	Inhibitor
Formula	$C_{20}H_{30}O_4$
MW	334.5
Purity	≥98% (TLC)
Solubility	Soluble in 100% ethanol, acetone or ethyl acetate.
Source	Synthetic.