# **KT5720**

#### **PKA** inhibitor

KT5720 is a hexylester derivative of K-252a and is a selective inhibitor of PKA (K  $_{\rm i}$  =0.056  $\mu M$  ). Cell permeable.

Citations: 12

View Online »

**Ordering Information** 

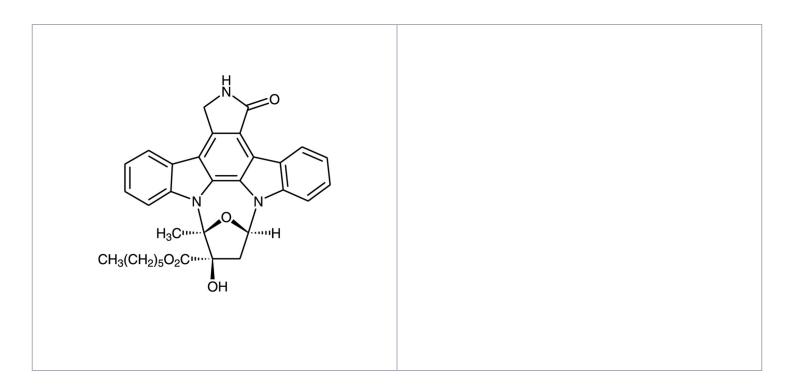
Order Online »

BML-EI199-0100

100µg

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. Stock solutions are stable for up

to 3 months at -20°C.

**Handling** Protect from light.

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

## Regulatory Status RUO - Research Use Only

### **Product Details**

**Appearance** White powder.

CAS 108068-98-0

Couple Target PKA

Couple Type Inhibitor

Formula  $C_{32}H_{31}N_3O_5$ 

MW 537.6

Purity ≥98% (HPLC)

**Solubility** Soluble in methanol: clear colorless solution at 5mg/ml.

Last modified: May 29, 2024

