# **Rac1** inhibitor

Cell-permeable, reversible, and selective Rac1 inhibitor.

NSC 23766 has been used to investigate the role of Rac1 in such diverse cellular functions as stem cell mobilization, epithelial cell migration, angiogenesis, leukemia cell migration and growth.

Citations: 3

View Online »

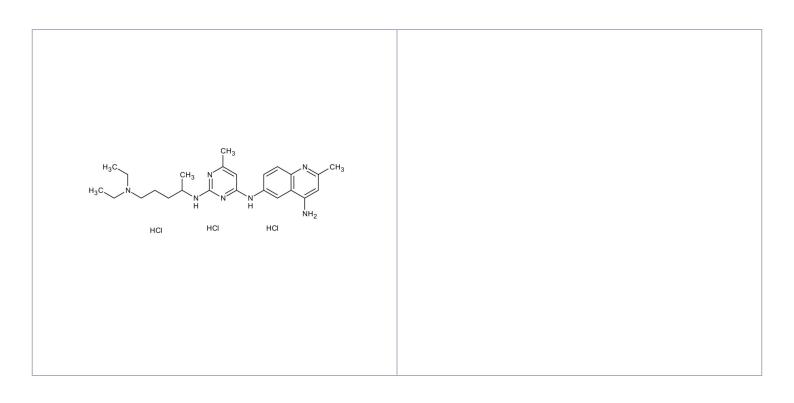
## **Ordering Information**

Order Online »

ENZ-CHM116-0005	5mg
ENZ-CHM116-0010	10mg

Manuals, SDS & CofA

View Online »



### **Handling & Storage**

**Use/Stability** As indicated on product label or CoA when stored as recommended.

Short Term Storage -20°C

**Long Term Storage** -80°C

**Shipping** Ambient Temperature

### Regulatory Status RUO - Research Use Only

#### **Product Details**

Alternative Name N6-[2-[[4-(Diethylamino)-1-methylbutyl]amino]-6-methyl-4-

pyrimidinyl]-2-methyl-4,6-quinolinediamine trihydrochloride

**Appearance** White powder solid.

**CAS** 733767-34-5

Couple Target Rac

Couple Type Inhibitor

Formula  ${\rm C}_{24}{\rm H}_{35}{\rm N}_7~.~3~{\rm HCI}$ 

**MW** 531

Purity ≥99% (HPLC)

**Solubility** Soluble in water.

Last modified: May 29, 2024

