# **SKF-525A. HCI**

#### **Cytochrome P450 inhibitor**

Cytochrome P-450 inhibitor. Inhibits cytochrome P-450-dependent arachidonate metabolism (90% at 50  $\mu$ M). Calcium antagonist. Blocks glibenclamide-sensitive K+ channels (IC $_{50}$ =4.4mM). Inhibits platelet thromboxane synthesis and stimulates endothelial cell prostacyclin production. Potentiates the effects of many drugs *in vivo*.

Citations: 13

View Online »

**Ordering Information** 

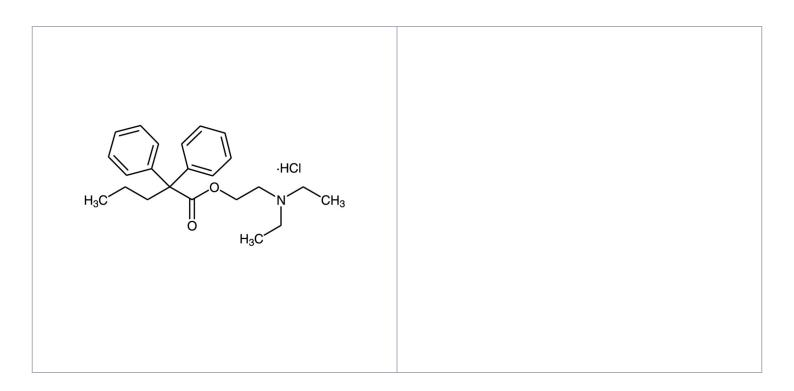
Order Online »

BML-EI105-1000

1g

Manuals, SDS & CofA

**View Online** »



### **Handling & Storage**

Use/Stability As indicated on product label or CoA when stored as recommended. Store, as supplied,

at room temperature for up to 1 year.

**Handling** Hygroscopic.

Long Term Storage Ambient

Shipping Ambient Temperature

## Regulatory Status RUO - Research Use Only

#### **Product Details**

Alternative Name Proadifen

**Appearance** White to off-white solid.

**CAS** 62-68-0

Couple Target Cytochrome P450, NOS, Potassium channel

Couple Type Blocker, Inhibitor

Formula  $C_{23}H_{31}NO_2$  . HCI

**Identity** Determined by 1H-NMR and MS.

**MW** 390.0

Purity ≥95% (TLC)

**Soluble** in water (25mg/ml) or 100% ethanol.

Last modified: May 29, 2024



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 info-

uk@enzolifesciences.com