

# SKF-525A . HCl

## 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<sup>+</sup> 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 »](#)

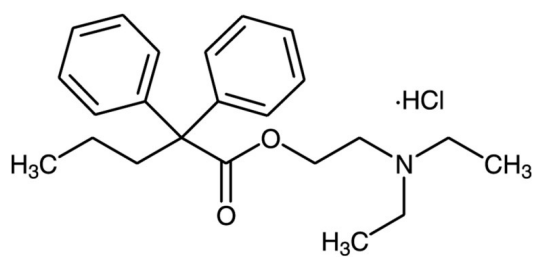
## Ordering Information

[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	$C_{23}H_{31}NO_2 \cdot HCl$
Identity	Determined by 1H-NMR and MS.
MW	390.0
Purity	≥95% (TLC)
Solubility	Soluble in water (25mg/ml) or 100% ethanol.



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)