

# Raloxifene . HCl

## Estrogen receptor inhibitor

A selective estrogen receptor modulator (SERM). Displays anti-estrogen action along with anti-androgen action. In clinical use for prevention of postmenopausal osteoporosis. Reduces circulating levels of IL-6 and TNF $\alpha$ . Inhibits prostate carcinogenesis in transgenic rat model. May be useful for the prevention of hormone-related cancers.

Citations: 2

[View Online »](#)

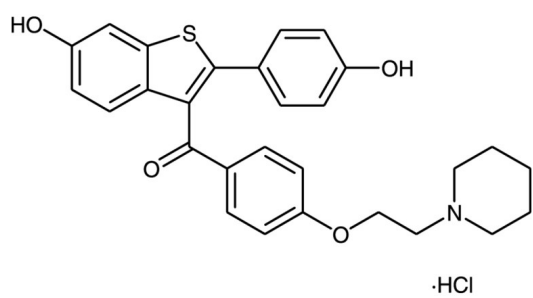
## Ordering Information

[Order Online »](#)

BML-GR243-0050	50mg
----------------	------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store reconstituted solutions at -20°C for up to three months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Appearance	Pale green solid.
CAS	82640-04-8
Couple Target	Estrogen receptor
Couple Type	Inhibitor, Ligand
Formula	$C_{28}H_{28}ClNO_4S$
Identity	Determined by $^1H$ -NMR and MS.
MW	510
Purity	≥98% (TLC)
Solubility	Soluble in methanol.



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)