

# A-803467

Na<sup>+</sup> channel blocker

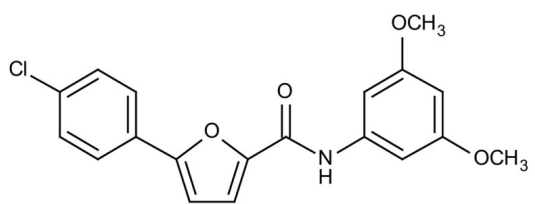
A-803467 is a potent and selective inhibitor of the TTX-resistant voltage-gated sodium channel Nav1.8 which is preferentially expressed on sensory neurons. IC<sub>50</sub> for human Nav1.8=8.0 nM (140 nM for rat). It attenuates neuropathic and inflammatory pain in various animal models.

Citations: 2

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Long Term Storage	Ambient
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Appearance	White solid.
CAS	944261-79-4
Couple Target	Sodium channel
Couple Type	Blocker
Formula	$C_{19}H_{16}ClNO_4$
Identity	Determined by NMR.
MW	357.8
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (15mg/ml) or 100% ethanol (2mg/ml).
Source	Synthetic.

Last modified: May 29, 2024



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-be@enzolifesciences.com](mailto:info-be@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)