

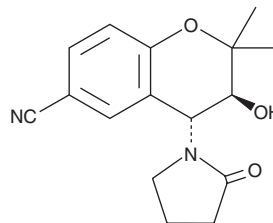
# PRODUCT INFORMATION



## (±)-Cromakalim

Item No. 14601

**CAS Registry No.:** 94470-67-4  
**Formal Name:** (3R,4S)-rel-3,4-dihydro-3-hydroxy-2,2-dimethyl-4-(2-oxo-1-pyrrolidinyl)-2H-1-benzopyran-6-carbonitrile  
**Synonym:** BRL 34915  
**MF:** C<sub>16</sub>H<sub>18</sub>N<sub>2</sub>O<sub>3</sub>  
**FW:** 286.3  
**Purity:** ≥98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

(±)-Cromakalim is supplied as a solid. A stock solution may be made by dissolving the (±)-cromakalim in the solvent of choice. (±)-Cromakalim is soluble in the organic solvent DMSO, which should be purged with an inert gas, at a concentration of approximately 50 mM.

### Description

(±)-Cromakalim is a prototypical activator of the ATP-sensitive potassium channel, SUR2-K<sub>ir</sub>6, that displays anti-ischemic (EC<sub>25</sub> = 8.9 μM) and cardioprotective (IC<sub>50</sub> = 0.03 μM) effects in rat hearts.<sup>1</sup> (±)-Cromakalim has also been shown to relax bladder smooth muscle *in vitro*.<sup>1</sup> It has been used to investigate potassium channel biology and as a scaffold to develop increasingly selective K<sup>+</sup> channel openers with clinical relevance.<sup>1</sup>

### Reference

1. Coghlan, M.J., Carroll, W.A., and Gopalakrishnan, M. Recent developments in the biology and medicinal chemistry of potassium channel modulators: Update from a decade of progress. *J. Med. Chem.* **44**(11), 1627-1653 (2001).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/13/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM