

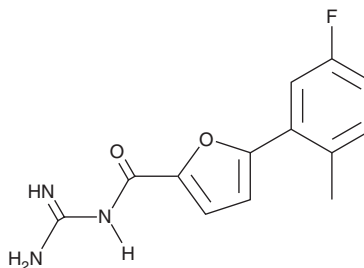
# PRODUCT INFORMATION



**KR-32568**

Item No. 23968

**CAS Registry No.:** 852146-73-7  
**Formal Name:** N-(aminoiminomethyl)-5-(5-fluoro-2-methylphenyl)-2-furancarboxamide  
**MF:** C<sub>13</sub>H<sub>12</sub>FN<sub>3</sub>O<sub>2</sub>  
**FW:** 261.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 316 nm  
**Supplied as:** A crystalline 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

KR-32568 is supplied as a crystalline solid. A stock solution may be made by dissolving the KR-32568 in the solvent of choice, which should be purged with an inert gas. KR-32568 is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

## Description

KR-32568 is an inhibitor of the sodium-hydrogen exchanger isoform-1 (NHE-1; IC<sub>50</sub> = 230 nM).<sup>1</sup> It restores cardiac contractile function *ex vivo* in an isolated ischemic rat heart model when used at a concentration of 10 μM. KR-32568 (0.3 mg/kg) reduces myocardial infarct size in a rat model of ischemia and reperfusion injury.

## Reference

1. Lee, S., Yi, K.Y., Hwang, S.K., *et al.* (5-Arylfuran-2-ylcarbonyl)guanidines as cardioprotectives through the inhibition of Na<sup>+</sup>/H<sup>+</sup> exchanger isoform-1. *J. Med. Chem.* **48**(8), 2882-2891 (2005).

### 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, 11/16/2022

## 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](http://WWW.CAYMANCHEM.COM)