

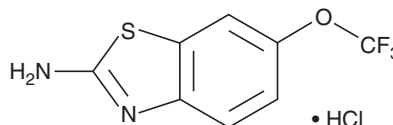
# PRODUCT INFORMATION



## Riluzole (hydrochloride)

Item No. 14577

**CAS Registry No.:** 850608-87-6  
**Formal Name:** 6-(trifluoromethoxy)-2-benzothiazolamine, monohydrochloride  
**Synonym:** PK 26124  
**MF:** C<sub>8</sub>H<sub>5</sub>F<sub>3</sub>N<sub>2</sub>OS • HCl  
**FW:** 270.7  
**Purity:** ≥99%  
**Supplied as:** A solid  
**Storage:** Room temperature  
**Stability:** ≥1 year



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

### Laboratory Procedures

Riluzole (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the riluzole (hydrochloride) in the solvent of choice. Riluzole (hydrochloride) is soluble in organic solvents such as DMSO, which should be purged with an inert gas. It is also soluble in water.

### Description

Riluzole is a benzothiazole derivative with anti-excitotoxic effects that acts by blocking the presynaptic release of glutamate, indirectly antagonizing glutamate receptors, and inactivating neuronal voltage-gated Na<sup>+</sup> channels (ED<sub>50</sub> = 2.3 μM).<sup>1</sup> Riluzole suppresses glutamate-induced seizures in rats at an ED<sub>50</sub> value of 3.2 mg/kg and displays neuroprotective effects in hypoxic animals at an ED<sub>50</sub> value of 4 mg/kg.<sup>1</sup> Formulations containing riluzole have been explored as therapeutics for slowing disease progression of amyotrophic lateral sclerosis.<sup>1</sup>

### Reference

1. Cheah, B.C., Vucic, S., Krishnan, A.V., *et al.* Riluzole, neuroprotection and amyotrophic lateral sclerosis. *Curr. Med. Chem.* **17(18)**, 1942-1959 (2010).

**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, 10/02/2017

### 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