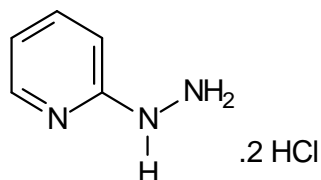


Product Data Sheet/Certificate of Analysis

2-Hydrazinopyridine · dihydrochloride

Storage: Store at room temperature



$C_5H_6N_3 \cdot 2HCl$ MW 182.5

Catalog Number:	S-2002	Unit Size:	
Lot Number:		Components:	2-Hydrazinopyridine.2HCl

Test	Specification	Result
Appearance (color)	Pale orange/brown	
Appearance (form)	Solid	
NMR:	>95%	

QC Release Date	Expiration Date:
Released By:	

Product Description

Quantitation of benzaldehyde moieties incorporated on proteins with 4FB.

Comments

Not for internal or external use in humans.

The products offered here are for research use only. Any commercial application will require a license from Solulink. The Solulink Conjugation System is patented and has multiple patents pending. Please contact Solulink for information regarding licensing information. Solulink products and methods may be covered by one or more of the following United States patents Nos. 6,686,461, 6,800,728, 7,102,024, 7,173,125, 7,462,689 and other pending patent applications. Information in this manual is subject to change without notice and does not constitute a commitment on the part of Solulink, Inc. It is supplied on an "as is" basis without any warranty of any kind, either explicit or implied. Information may be changed or updated in this manual at any time. This document may not be copied, transferred, reproduced, disclosed, or duplicated, in whole or in part, without the prior written consent of Solulink, Inc. This documentation is proprietary information and protected by the copyright laws of the United States and international treaties. The manufacturer of this documentation is Solulink, Inc.

