

# Propidium Iodide

Product number: CSK-0118A

Lot Number: 3408418

Test	Specification	Result
Absorption Spectrum $\lambda$ (potassium phosphate buffer)(nm) $\epsilon$ (493)	$493 \pm 2$	Passed
	$5,900 \pm 500$	Passed
HPLC	$\geq 98\%$	Passed
TLC	$\geq 95\%$	Passed

Manufacturing date: 05-May-2025

Document version: B1

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.