



BIOSCIENCES®

G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: N-Tosyl-L-alanine 3-indoxyl ester**Catalog No:** RC-796**Release Date:** 8/17/20**Size:** 1g**Retest Date:** 6/2023**Lot No:** 201708**Storage:** -20°C

Specifications

CAS Number: 75062-54-3**Appearance:** White to off-white crystalline powder

Test	Specification	Result
Color of solution (1% in methanol)	Colorless	Colorless
Turbidity (1% in methanol)	≤4.0 NTU	0.4 NTU
Water Content (K. Fischer)	≤1.0%	0.3%
Melting Point	75-79°C	78°C
Rotation	-57.0 to -53.0	-56.3
Purity (HPLC)	≥98.0 area-%	99.8% area-%
TLC	Corresponds	PASS
FT-IR Spectrum	Corresponds	PASS

This document has been produced electronically and is valid without signature

Date of Print: September 14, 2020**G-Biosciences/Geno Technology, Inc.**

Rev 10/25/18 BP

**think proteins! think G-Biosciences www.GBiosciences.com**