



## Certificate of Analysis / Data Sheet

**Product:** Aprotinin

**Catalog No:** 786-1245

**Size:** 25mg

**Lot No:** 203101

**Release Date:** 1/31/20

**Retest Date:** 1/31/22

**Storage:** 4°C

### Specifications

**CAS Number:** 9087-70-1

**Appearance:** White to off-white hygroscopic powder

Test	Specification	Result
Solubility	Clear, colorless (2mg/ml water)	PASS
Purity (HPLC)	≥ 95.0%	99.5%
Activity	3.0-8.0 TIU/mg solid	4.9 TIU/mg solid
Protein content	≥ 0.9 mg protein/mg solid	0.91mg protein/mg solid
pH	5.0-7.0 (5mg/ml water)	6.1
Loss on Drying	≤ 6.0%	5.2%

This document has been produced electronically and is valid without signature

**Date of Print:** February 5, 2020

**G-Biosciences/Geno Technology, Inc.**

Rev 10/25/18 BP

