



**Data Sheet /**  
**Certificate of Analysis**

**Product:** Casein-Resorufin Protease Substrate

**Catalog #:** 786-321

**Size:** 5mg

**Lot #:** 100206

**Description:** Resorufin-casein Protease Substrate is a milk protein (casein) labeled with resorufin [N-(resorufin-4-carbonyl) piperidine-4-carbonic acid]. It is a general protease substrate and can be used for sensitive detection of protease contamination in the test sample, in both colorimetric and fluorometric methods. The treatment of the substrate resorufin-casein with protease(s) results in the release of resorufin-labeled peptides, which are not precipitated by trichloroacetic acid. The concentration of the labeled peptides in the supernatant is proportional to the proteolytic activity present in the sample.

**Tests:**

**Appearance:** Lyophilized

**Form:** Solid

**Solubility:** 4mg/ml in DI water

**Color:** Purple red

**Storage:** Store at -20<sup>0</sup>C. After its reconstitution, aliquots of smaller volume can be stored at 4<sup>0</sup>C for short term storage (2-3 days) and at -20<sup>0</sup>C for long term storage (2-3 months).

This document has been produced electronically and is valid without signature

**Date of Print:** December 1, 2010

**G-Biosciences / Genotech**

92 Weldon Parkway, St. Louis, MO 63043, USA

Fax: 314-991-1504

Casein-Resorufin Protease Substrate CofA/09.14.09/IA/Common3



**think proteins! think G-Biosciences!**

**[www.GBiosciences.com](http://www.GBiosciences.com)**