

Certificate of Analysis

EarlyTox™ Caspase-3/7 R100 Assay Explorer Kit

R8346

Lot #	1757737
--------------	----------------

R8346

	Specification	This Lot
Cell-based Caspase-3 Assay Test	Meets specification	pass

Expiry Date	Up to 6 months from receipt when stored at -20°C
--------------------	---

This product has been manufactured to ISO9001:2008 Quality Assurance guidelines

For research use only Store at -20°C

Signed: *Erwin Ramirez* Date: 24-May-2016

Production Chemist