

Certificate of Analysis

EarlyTox™ Caspase-3/7 NucView™ 488 Assay Explorer Kit R8350

Lot #	1757736
--------------	---------

R8350A, NucView™ 488 Caspase-3 Substrate, 1mM in PBS

	Specification	This Lot
Max. Abs. (MeOH) (nm)	503±3	501
TLC	≥95%	95%
OD at λmax (DMSO)	63±3	63.2

R8348B, Masking Reagent

	Specification	This Lot
Moisture Test (weight loss)	<3%	2.67%

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