

# Certificate of Analysis

## LiveDetect Green

### K8300

<b>Batch #</b>	<b>LDG1212A</b>
----------------	-----------------

<b>Full Description</b>	Fluorescent cytoplasmic dye for the detection of viable colonies of mammalian cells in semi solid culture.
<b>Components</b>	No material of animal origin is used in the manufacture of this product.
<b>Appearance</b>	Clear solution
<b>Concentration</b>	1mM
<b>Format</b>	Liquid (frozen for storage)
<b>Volume</b>	50µl in 5ml spray bottle
<b>Functionality Test</b>	Detection of green fluorescence (FITC channel) of viable mammalian colonies in semi solid media after >18 hours incubation.
<b>Cell Line Used</b>	Hybridoma Cell Line SJK-287-38 ECACC Cat # 90112228
<b>Solvent</b>	DMSO (Sigma Product: D2650, Lot: RNBB8135)
<b>Sterility</b>	No bacterial or fungal growth after 72 hours incubation
<b>Preservatives/Stabilisers</b>	None
<b>Date of manufacture</b>	13-Dec-2012
<b>Expiry Date</b>	13-Jun-2013

This product has been manufactured to ISO9001:2008 certification

**For research use only    Store at -20°C**

Signed:

Date: 13-Dec-2012

*Christopher Lee*

**Production Chemist**