

# Certificate of Analysis

## Calibeads

### K8150

<b>Batch #</b>	<b>CBA0714A</b>
----------------	-----------------

<b>Full Description</b>	Uniform fluorescent polystyrene microspheres in methylcellulose. Molecular Devices Calibeads are microspheres that fluoresce at four distinct wavelengths corresponding to the CFP, FITC/GFP, Rhodamine/PE and Cy5 filter sets of ClonePix FL / ClonePix2
<b>Appearance</b>	Clear viscous liquid
<b>Components</b>	200µm polystyrene spheres
<b>Volume</b>	100ml in semi solid media
<b>Concentration</b>	ca. 10-25 beads/ml
<b>Sterility</b>	No bacterial or fungal growth after 72 hours incubation
<b>Preservatives/Stabilisers</b>	None
<b>Buffer</b>	1 x PBS pH 7.2
<b>Functionality Test</b>	Tested and validated in functional assay on ClonePix FL / ClonePix2
<b>Date of manufacture</b>	27-Jul-2014
<b>Expiry Date</b>	Up to 3 years from receipt when stored in the dark at 2°C – 8°C

This product has been manufactured to ISO9001:2008 certification

**For research use only    Store at 2°C – 8°C    Light sensitive**

Signed:

Date: 31-Jul-2014



**Production Biologist**