

Certificate of Analysis

Calibeads

K8150

Batch #	CBA0219A
----------------	-----------------

Full Description	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
Appearance	Clear viscous liquid
Components	200µm polystyrene spheres
Volume	100ml in semi solid media
Concentration	ca. 10-25 beads/ml
Sterility	No bacterial or fungal growth after 72 hours incubation
Preservatives/Stabilisers	None
Buffer	1 x PBS pH 7.2
Functionality Test	Tested and validated in functional assay on ClonePix FL / ClonePix2
Expiry Date	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:

Kieu-Khiem Nguyen

Date: 04-FEB-2019

Production Biologist