



## CERTIFICATE OF ANALYSIS

Product Name: Clone Select™ Imager Reference Plate

Part Number: 5030520

Lot #: CSR0515A

Storage: Store at Room Temp.

Expiry Date: Up to 12 months from receipt

Overall Result:	<b>PASS</b>
This document certifies that the product defined below has achieved a "PASS" on all required performance, functional and/or safety verification tests established by Molecular Devices in accordance with applicable regulatory requirements as summarized below.	

### Results

		Specification
<b>CELL CONFLUENCE</b>	<b>PASS</b>	Highest concentration wells (Column 1 & 2) Confluence of >70%. Lowest concentration wells (Column 11 & 12) Confluence of <10%.
<b>CELL NUMBER</b>	<b>PASS</b>	Cell number within tolerance
<b>CloneSelect Imager</b> Inspection	<b>PASS</b>	Cells uniformly distributed
<b>MISCELLANEOUS INFORMATION</b>	<b>PASS</b>	Material lot Number

Approved by: \_\_\_\_\_ *Ervin Ramirez* \_\_\_\_\_

Date: \_\_\_\_\_ 22-May-2015 \_\_\_\_\_

Production Biologist