

Product Name: Clone Select™ Imager Reference Plate

Part Number: 5030520

Lot #: 073014

Storage: Store at 2 - 8 °C

Overall Result:	<b>PASS</b>
<p>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.</p>	

## 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 of
<b>CloneSelect Imager Inspection</b>	<b>PASS</b>	Cells uniformly distributed
<b>MISCELLANEOUS INFORMATION</b> Material lot Number	<b>PASS</b>	

Approved by: *Christopher Lee* 30-Jul-2014