

Product Name: Clone Select™ Imager Reference Plate

Part Number: 5030520

Lot #: CSR0914A

Storage: Store at 2 - 8 °C

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

Approved by: Christopher Lee 11-Sep-2014