

Certificate of Analysis

Product name OncoSpot™ Wild type HCT116 Genomic DNA Reference Standard
Product Description Wild type HCT116 Genomic DNA Reference Standard
Cat. No. RM001
Lot. No. 325EQ0101
Unit Size 5ug (50ng/ul)

TEST	TEST METHOD	ACCEPTANCE CRITERIA	PASS/FAIL
Genotype	Sanger sequencing of locus specific PCR	Correct base alteration at chromatograph matching the gDNA zygosity	PASS
Quality	Agarose gel electrophoresis	Bright band of high molecular weight	PASS
Quantification	Qubit 3.0 Fluorometer	50±5 ng/μL	PASS

Conclusion This lot passes the Quality Control specifications.

Signed by: 
Quality Inspector

GeneCopoeia, Inc.

Address: 9620 Medical Center Drive, #101
Rockville, MD 20850
USA

Phone: 301-762-0888 **Fax:** 301-762-3888

Website: www.genecopoeia.com

E-Mail: inquiry@genecopoeia.com; support@genecopoeia.com