



CGH Labeling Kit (ENZ-42671): Cited Samples

Cells

SAMPLE TYPE	SAMPLE ORIGIN	TYPE OF ARRAY	PLATFORM	REFERENCES
Cells	Amniotic fluid	4x44K Agilent array CGH	Agilent Technologies	A.L. Mosca-Boidron, et al. (2013)
Cells	Blood	180K custom Oxford array CGH	Oxford Gene Technology	J. Lundin, et al. (2019)
Cells	Blood	60K Agilent array CGH	Agilent Technologies	E. Yokoyama, et al. (2018)
Cells	Blood	180K custom Oxford array CGH	Oxford Gene Technology	S. Arkani, et al. (2018)
Cells	Blood	Agilent array CGH	Agilent Technologies	D.N. Senanayake, et al. (2015)
Cells	Blood	BAC array CGH	Agilent Technologies	S. Struski, et al. (2007)
Cells	Blood	4x44K Agilent array CGH	Agilent Technologies	B. Grisart, et al. (2008)
Cells	Blood	4x44K Agilent array CGH	Agilent Technologies	J.W. Ahn, et al. (2010)
Cells	Blood	105K Agilent array CGH	Agilent Technologies	A.B.P. van Kuilenburg, et al. (2010)
Cells	Blood	4x44K Agilent array CGH	Agilent Technologies	A. Hills, et al. (2010)
Cells	Blood	4x44K Agilent array CGH	Agilent Technologies	S. Hoang, et al. (2011)
Cells	Blood	4x180K Agilent array CGH	Agilent Technologies	F.S. van Dijk, et al. (2010)
Cells	Blood	4x180K Agilent array CGH	Agilent Technologies	J.B. van de Meerakker, et al. (2011)
Cells	Blood	2x105K Oxford array CGH	Agilent Technologies	D.Q. Barge-Schaapveld, et al. (2011)
Cells	Blood	4x180K Agilent array CGH	Agilent Technologies	K. Szakson, et al. (2012)
Cells	Blood	Exon-level array CGH	Agilent Technologies	S. Aradhya, et al. (2012)
Cells	Blood	4x180K Agilent array CGH	Agilent Technologies	E. Markljung, et al. (2012)
Cells	Blood	4x180K Agilent array CGH	Agilent Technologies	J. Aury-Landas, et al. (2013)
Cells	Blood	60K Agilent array CGH	Agilent Technologies	M. Tropeano, et al. (2013)
Cells	Blood	1M Oxford array CGH	Agilent Technologies	A. Norling, et al. (2013)
Cells	Blood	180K Oxford array CGH	Agilent Technologies	J. Winberg, et al. (2014)
Cells	Blood	60K Agilent array CGH	Agilent Technologies	J.W. Ahn, et al. (2014)
Cells	Blood	180K and 244K Agilent array CGH	Agilent Technologies	A. Adegbola, et al. (2015)
Cells	Blood	4x180K Agilent array CGH	Agilent Technologies	J.P. Vieira, et al. (2015)
Cells	Gastric cancer cell lines	4x44K Agilent array CGH	Agilent Technologies	T.E. Buffart, et al. (2009)
Cells	Huntington's disease human embryonic stem cells	4x44K Agilent array CGH	Agilent Technologies	L. Jacquet, et al. (2015)
Cells	Induced pluripotent stem cells	60K Agilent array CGH	Agilent Technologies	T. Yagi, et al. (2011)
Cells	Lymphoblastoid cell lines	180K and 244K Agilent array CGH	Agilent Technologies	A. Adegbola, et al. (2015)
Cells	Neuroblastoma cell lines	BAC array CGH	Molecular Devices	H. Auer, et al. (2007)
Cells	Induced pluripotent stem cell lines	4x44K Agilent array CGH	Agilent Technologies	A. Petrova, et al. (2016)
Cells	Sperm	Large-insert clone whole-genome tilepath array	Agilent Technologies	D.J. Turner, et al. (2008)

Complete citation can be found on product's webpage

Global Headquarters
ENZO LIFE SCIENCES, INC.
 Ph: 800.942.0430
 info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES (ELS) AG
 Ph: +41 61 926 8989
 info-eu@enzolifesciences.com

Belgium, The Netherlands & Luxembourg
 Ph: +32 3 466 0420
 info-be@enzolifesciences.com

France
 Ph: +33 472 440 655
 info-fr@enzolifesciences.com

Germany
 Ph: +49 7621 5500 526
 info-de@enzolifesciences.com

UK & Ireland
 Ph: 0845 601 1488 (UK customers)
 Ph: +44 1392 825900
 info-uk@enzolifesciences.com

For local distributors and detailed product information visit us online: www.enzolifesciences.com

For Research Only. Not for Use in Diagnostic Procedures.

FFPE Tissue

SAMPLE TYPE	SAMPLE ORIGIN	TYPE OF ARRAY	PLATFORM	REFERENCES
FFPE tissue	Adenoid cystic carcinoma	4x180K Agilent array CGH	Agilent Technologies	A.F. Costa, et al. (2014)
FFPE tissue	Breast cancer tissue	135K NimbleGen array CGH	Agilent Technologies	P.C. Schouten, et al. (2015)
FFPE tissue	Diffuse large B-cell lymphoma	Custom Agilent array CGH	Agilent Technologies	L.M. Dias, et al. (2015)
FFPE tissue	Epithelial ovarian tumors	4x180K Agilent array CGH	Agilent Technologies	M.M. Kamieniak, et al. (2013)
FFPE tissue	Melanoma and melanocytic spitz nevi	BAC array CGH	Molecular Devices	L. Ali, et al. (2010)
FFPE tissue	Microdissected intraductal and invasive breast cancer	BAC array CGH	Molecular Devices	S. Liao, et al. (2012)
FFPE tissue	Polypoid and flat colorectal adenoma	4x44K Agilent array CGH	Agilent Technologies	Q.J. Voorham, et al. (2012)
FFPE tissue	Primary ovarian tumor tissue	2x105K Oxford array CGH	Agilent Technologies	A. Zaal, et al. (2012)
FFPE tissue	Stage II colon cancer samples	4x44K Agilent array CGH	Agilent Technologies	R.P. Brosens, et al. (2011)

Other Samples

SAMPLE TYPE	SAMPLE ORIGIN	TYPE OF ARRAY	PLATFORM	REFERENCES
ChIP DNA	ChIP-seq DNA library	60K Agilent array CGH	Agilent Technologies	D. Schmidt, et al. (2008)
ChIP DNA	Drosophila S2 cells	D.melanogaster NimbleGen array CGH	NimbleGen	H. Mason-Suares, et al. (2013)
Fresh tissue	Mouse tail	BAC array CGH	Molecular Devices	Z. Li, et al. (2007)
Frozen cells	P.gingivalis	P.gingivalis strain W83 array CGH	Agilent Technologies	J. Brunner, et al. (2010)
Frozen tissue	Primary uveal Melanoma	BAC array CGH	Molecular Devices	L.A. Worley, et al. (2007)
Frozen tissue	Heart, liver, lung and spleen tissue	180K Oxford array CGH	Agilent Technologies	J. Winberg, et al. (2014)

For more information, go to:

www.enzolifesciences.com/CGH