

OncoScan™ CNV Assay for Research

Catalog Numbers 902695

Pub. No. MAN0027555 Rev. D00



WARNING! Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from thermofisher.com/support.

Product use

The Applied Biosystems™ OncoScan™ CNV Assay for Research includes Applied Biosystems™ OncoScan™ CNV Arrays and the Applied Biosystems™ OncoScan™ CNV Reagent Kit for Research. The OncoScan™ CNV Assay for Research detects genome-wide copy number gains and losses and loss of heterozygosity (LOH), including copy neutral LOH, using limited DNA from Formalin-Fixed Paraffin-Embedded (FFPE) samples of various ages.

See thermofisher.com for a list of supporting manuals for assay and data analysis procedures.

Instructions for use

See the *OncoScan™ CNV Assay User Guide* (Pub. No. [MAN0027927](http://thermofisher.com)) and the *OncoScan™ CNV Assay—7 Samples Quick Reference* (Pub. No. [MAN0027190](http://thermofisher.com)) for procedures on DNA sample preparation, sample hybridization, fluidics setup, array scanning, and data analysis instructions.

OncoScan™ CNV Reagent Kit for Research components (Cat. No. [902692TS](http://thermofisher.com))

Module	Part No. ^[1]
OncoScan™ CNV Buffer C <ul style="list-style-type: none"> • Buffer C 	902687 <ul style="list-style-type: none"> • 902696
OncoScan™ CNV Copy Number Probe Mix 1.0 and Controls <ul style="list-style-type: none"> • Positive Control (12 ng/μL) • Negative Control • Copy Number Probe Mix 1.0 • Buffer A 	902688 <ul style="list-style-type: none"> • 902249 • 902250 • 902248 • 902246
OncoScan™ CNV Gap Fill and 1st Stage PCR <ul style="list-style-type: none"> • Buffer A • Gap Fill Enzyme Mix • SAP, Recombinant (1 U/μL) • dNTP Mix (A/T) • dNTP Mix (G/C) • Nuclease-Free Water • Exo Mix • Cleavage Buffer • Cleavage Enzyme (2 U/μL) • PCR Mix • Taq Polymerase (5 U/μL) 	902689 <ul style="list-style-type: none"> • 902246 • 902252 • 902251 • 902254 • 902255 • 902253 • 902256 • 902257 • 902258 • 902259 • 902260
OncoScan™ CNV 2nd Stage PCR and Post PCR Processing <ul style="list-style-type: none"> • PCR Mix • Taq Polymerase (5 U/μL) • Buffer B • HaeIII Enzyme (10 U/μL) • Exo I Enzyme (20 U/μL) • Nuclease-Free Water • Hybridization Mix 	902690 <ul style="list-style-type: none"> • 902259 • 902260 • 902261 • 902262 • 902263 • 902253 • 902264
OncoScan™ CNV Stain Reagents <ul style="list-style-type: none"> • Stain 1 • Stain 2 • Array Holding Buffer 	902691 <ul style="list-style-type: none"> • 902265 • 902266 • 901733
<ul style="list-style-type: none"> • Wash A (Individual bottle) • Wash B (Individual bottle) 	<ul style="list-style-type: none"> • 901680 • 901681

^[1] Part numbers are for identification purposes only and cannot be ordered separately.

Storage, handling, and stability

The reagents that are contained in the OncoScan™ CNV Assay for Research are shipped at different temperatures and must be stored at the recommended conditions.

The following modules must be stored at -20°C:

- OncoScan™ CNV Buffer C
- OncoScan™ CNV Copy Number Probe Mix 1.0 & Controls
- OncoScan™ CNV Gap Fill and 1st Stage PCR
- OncoScan™ CNV 2nd Stage PCR

The following modules must be stored at 2–8°C:

- OncoScan™ CNV Stain Reagents
- OncoScan™ CNV Arrays

The following individual bottles must be stored at room temperature:

- Wash A
- Wash B

See the expiration date on the package label. Do not use probe arrays or reagents components after the expiration date.

IMPORTANT! Store the unopened kits in rooms appropriate to the workflow. The reagent subkits are separated into modules to support the recommended workflow and minimize the chance of contamination during use.

Ordering information

Cat. No.	Product name	Description
902695	OncoScan™ CNV Assay for Research, consists of: <ul style="list-style-type: none">• OncoScan™ CNV Reagent Kit for Research• 48 OncoScan™ CNV Arrays	Sufficient for 24 samples
902693	OncoScan™ CNV Assay Training Kit, consists of: <ul style="list-style-type: none">• OncoScan™ CNV Reagent Kit for Research• 36 OncoScan™ CNV Arrays• Training controls	Sufficient for 18 samples
902692TS	OncoScan™ CNV Reagent Kit for Research	Sufficient for 24 samples
902694	OncoScan™ CNV Arrays	48 arrays

Limited product warranty

Life Technologies Corporation and its affiliates warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have questions, contact Life Technologies at www.thermofisher.com/support.



Thermo Fisher Scientific Baltics UAB | V.A. Graiciuno 8, LT-02241 | Vilnius, Lithuania

Revision history: Pub. No. MAN0027555 D00

Revision	Date	Description
D00	10 April 2024	<ul style="list-style-type: none">• Initial release in Thermo Fisher Scientific document control system. Supersedes legacy Affymetrix publication number 703329.• Added "for Research" to product name and title.
3	17 August 2017	Remove the word "Manufacturer" next to symbol on page 2.
2	1 August 2017	Updated product name. Updated document to current template.
1	31 March 2015	Baseline for revision history.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2015-2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.