

OncoScan™ CNV Plus Assay for Research

Catalog Numbers 902293

Pub. No. MAN0027540 Rev. E00



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](https://www.thermofisher.com/support).

Product use

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

OncoScan™ CNV Plus Assay for Research—same copy number coverage as the OncoScan™ CNV Assay for Research plus a somatic mutation panel covering 74 mutations in nine genes (BRAF, KRAS, EGFR, IDH1, IDH2, PTEN, PIK3CA, NRAS, TP53).

See [thermofisher.com](https://www.thermofisher.com) for a list of supporting manuals for assay and data analysis procedures.

Instructions for use

See the *OncoScan™ CNV Plus Assay User Guide* (Pub. No. [MAN0027850](https://www.thermofisher.com)) and the *OncoScan™ CNV Plus Assay—7 Samples Quick Reference* (Pub. No. [MAN0029389](https://www.thermofisher.com)) for procedures on DNA sample preparation, sample hybridization, fluidics setup, array scanning, and data analysis instructions.

OncoScan™ CNV Plus Reagent Kit for Research components (Cat. No. [902294](https://www.thermofisher.com))

Module	Part No. ^[1]
OncoScan™ CNV Plus Somatic Mutation Probe Mix 1.0 <ul style="list-style-type: none"> OncoScan™ CNV Plus Somatic Mutation Probe Mix 1.0 	902272 <ul style="list-style-type: none"> 902247
OncoScan™ CNV Plus Somatic Mutation Probe Mix 1.0 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 	902268 <ul style="list-style-type: none"> 902249 902250 902248 902246
OncoScan™ CNV Plus 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) 	902269 <ul style="list-style-type: none"> 902246 902252 902251 902254 902255 902253 902256 902257 902258 902259 902260
OncoScan™ CNV Plus 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 	902270 <ul style="list-style-type: none"> 902259 902260 902261 902262 902263 902253 902264
OncoScan™ CNV Plus Stain Reagents <ul style="list-style-type: none"> Stain 1 Stain 2 Array Holding Buffer 	902271 <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] For identification purposes only and cannot be ordered separately.

Storage, handling, and stability

The reagents that are contained in the OncoScan™ CNV Plus 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 Plus Somatic Mutation Probe Mix 1.0
- OncoScan™ CNV Plus Copy Number Probe Mix 1.0 and Controls
- OncoScan™ CNV Plus Gap Fill and 1st Stage PCR
- OncoScan™ CNV Plus 2nd Stage PCR and Post PCR Processing

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

- OncoScan™ CNV Plus Stain Reagents
- OncoScan™ CNV Plus 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
902293	OncoScan™ CNV Plus Assay for Research, consists of: <ul style="list-style-type: none">• OncoScan™ CNV Plus Reagent Kit for Research• 48 OncoScan™ CNV Plus Arrays	Sufficient for 24 samples
902305	OncoScan™ CNV Plus Assay Training Kit, consists of: <ul style="list-style-type: none">• OncoScan™ CNV Plus Reagent Kit for Research• 36 OncoScan™ CNV Plus Arrays• Training controls	Sufficient for 18 samples
902294	OncoScan™ CNV Plus Reagent Kit for Research	Sufficient for 24 samples
902292	OncoScan™ CNV Plus Array	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. MAN0027540 E00

Revision	Date	Description
E00	17 April 2024	<ul style="list-style-type: none">• Initial release in Thermo Fisher Scientific document control system. Supersedes legacy Affymetrix publication number 703178.• Added "for Research" to product name and title.
4	17 August 2017	Remove the word "Manufacturer" next to symbol on page 2.
3	1 August 2017	Updated document to current template.
2	11 July 2013	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.

©2013-2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.