

**For In Vitro Diagnostic Use**

HD832-IVD Mimix™ OncoSpan FFPE Reference Standard

**Intended Use**

The Mimix™ OncoSpan FFPE reference standard is a commutable control material comprising a formalin-fixed paraffin-embedded (FFPE) curl derived from human cancer cell lines, containing more than 380 variants across 152 cancer genes. It is intended for qualitative and/or quantitative monitoring of next-generation sequencing (NGS) or droplet digital polymerase chain reaction (ddPCR) assays designed to detect somatic mutations in genomic DNA (gDNA) from human samples for *in vitro* diagnostic use.

The gDNA obtained from OncoSpan FFPE Reference Standard can be used in NGS or ddPCR workflow, test performance, assay variation and helps identify increases in random or systematic errors. This product is for professional laboratory use only.

**Summary and Explanation**

NGS assays are widely adopted by researchers as well as clinical and diagnostic laboratories to analyze the genetic make-up of patient-derived samples, making it useful to gain insightful data on common or rare diseases<sup>1</sup>. Providing high-throughput data, NGS helps identify genetic differences in patients or any genetic changes introduced as certain diseases progress<sup>2</sup>. Similarly, digital PCR workflows, including ddPCR, are commonly used to enhance traditional PCR performance by allowing a significantly increased number of reactions to be performed and analyzed real time<sup>3</sup>.

The Mimix™ OncoSpan FFPE Reference Standard enables direct analysis of patient-derived DNA samples against a wide range of cancer-related somatic mutations for fast and accurate results as part of diagnostic tests and scientific research. The product is designed to fit into NGS and ddPCR workflows to help accurately analyze patient-derived samples.

OncoSpan FFPE standard is a curl mimicking the patient tissue sections, sectioned from an FFPE block derived from well-established cell lines via cell pellet collection and blending followed by fixation in 4% paraformaldehyde (PFA) and paraffin embedding. Each FFPE block contains

approximately  $3 \times 10^8$  cells, which corresponds to approximately  $3.5 \times 10^5$  cells per section. The standard contains more than 380 variants across 152 key cancer genes, with variant allele frequencies ranging from 1% to 92.5% with 52 variants present at  $\leq 20\%$  allelic frequency for limit of detection determination of your assay. This diversity allows for comprehensive assessment of assay performance and quantitative accuracy.

The Mimix™ OncoSpan FFPE Reference Standard is a multiplex standard covering the mutations on the following genes: ABL1, AKT1, AKT2, ALK, APC, AR, ARID1A, ATR, ATRX, AXL, BARD1, BCL6, BLM, BMPR1A, BRAF, BRCA1, BRCA2, BTK, BUB1B, CARD11, CCND1, CCND3, CCNE1, CD79B, CDH2, CDK12, CDK4, CEP57, CFH, CREBBP, CSF1R, CTNNB1, DDR2, DIS3L2, DNMT3A, EGFR, EML4, EP300, EPCAM, ERBB2, ERBB3, ERCC1, ERCC2, ERCC4, ERCC5, ERG, ETS1, ETV4, EWSR1, EXT1, FANCA, FANCD2, FANCE, FANCG, FANCI, FANCM, FBXW7, FGF10, FGF2, FGF3, FGF6, FGFR1, FGFR3, FLCN, FLI1, FLT1, FLT3, FZR1, GATA2, GATA3, GEN1, GNA11, GNAS, HNF1A, HRAS, IDH2, IDH4, JAK1, JAK2, JAK3, KDR, KIT, KRAS, LDLR, MAGI1, MAP2K1, MAP2K2, MAX, MDM4, MED12, MET, MLH2, MLLT3, MMAB, MRE11, MSH4, MSH6, MTOR, NBN, NF1, NFE2L2, NOTCH1, NOTCH4, NOTCH6, NRAS, NRG1, NTRK1, NTRK3, PDGFRA, PDGFRB, PIK3CA, PIK3CD, PIK3CG, PIK3R1, PMS2, PPARG, PPP2R2A, PRKAR1A, PROC, PTCH2, PTPN11, RAD51B, RAD54L, RAF1, RB1, RBM45, RECQL4, RET, RHBDF2, ROS1, RPS6KB1, SDHB, SF3B1, SF3B2, SLTM, SLX4, SMARCB1, SMO, SMOX, STK11, TERT, TET2, TFRC, TP53, TP53BP1, TSC1, TSC2, WRN, XPA, XPC, ZNF395.

**Principles of Operation**

The Mimix™ OncoSpan FFPE Reference Standard contains a single curl suitable for gDNA extraction, derived directly from well-characterized cell lines to mimic patient samples. No synthetic DNA is incorporated into this product. The FFPE curl is intended to be integrated into the customer's workflow for DNA extraction. Every batch of OncoSpan FFPE has 25 variants confirmed by ddPCR (Table 1), the other 355 variants are covered by exome sequencing. This provides an accurate and reliable set for comparison to any NGS assay's performance.

In addition to ddPCR and NGS, further quality control of this product was performed via agarose gel electrophoresis and fluorometric analysis.

Table 1: Verified variants and the corresponding allelic frequencies (AF) confirmed by ddPCR.

| Chromosome       | Gene   | Variant              | Expected AF | Acceptance Range |
|------------------|--------|----------------------|-------------|------------------|
| chr2 (29416025)  | ALK    | N/A (Ins)            | 10.0%       | 8.00 – 12.00%    |
| chr5 (112175770) | APC    | p.T1493T             | 35.0%       | 31.50 – 38.50%   |
| chr7 (140453136) | BRAF   | p.V600E              | 10.70%      | 7.49 – 13.91%    |
| chr13 (32913558) | BRCA2  | p.K1691fs*15         | 32.5%       | 29.25 – 35.75%   |
| chr3 (41266101)  | CTNNB1 | p.S33Y               | 32.5%       | 29.25 – 35.75%   |
| chr3 (41266133)  | CTNNB1 | p.S45del             | 10.0%       | 8.00 – 12.00%    |
| chr7 (55241707)  | EGFR   | p.G719S              | 24.5%       | 22.10 – 27.00%   |
| chr7 (55249071)  | EGFR   | p.T790M              | 1.0%        | 0.60 – 1.20%     |
| chr7 (55259515)  | EGFR   | p.L858R              | 3.0%        | 1.68 – 3.92%     |
| chr7 (55242464)  | EGFR   | p.E746_A750 delELREA | 2.0%        | 1.30 – 2.50%     |
| chr7 (55249063)  | EGFR   | p.Q787Q              | 15.0%       | 12.00 – 18.00%   |
| chr4 (153244155) | FBXW7  | p.S668fs*39          | 32.5%       | 29.25 – 35.75%   |
| chr13 (28578214) | FLT3   | p.P986fs*>8          | 10.0%       | 8.00 – 12.00%    |
| chr4 (55599321)  | KIT    | p.D816V              | 10.0%       | 8.00 – 12.00%    |
| chr4 (55602765)  | KIT    | p.L862L              | 7.5%        | 6.00 – 9.00%     |
| chr12 (25398281) | KRAS   | p.G13D               | 15.0%       | 12.00 – 18.00%   |
| chr12 (25398284) | KRAS   | p.G12D               | 6.0%        | 5.00 – 7.60%     |
| chr7 (116436022) | MET    | p.A1357A             | 7.0%        | 5.60 – 8.40%     |
| chr7 (116339847) | MET    | p.L238fs*25          | 7.0%        | 5.60 – 8.40%     |
| chr9 (139409754) | NOTCh1 | p.P668S              | 30.0%       | 27.00 – 33.00%   |
| chr1 (115256530) | NRAS   | p.Q61K               | 12.5%       | 10.00 – 15.00%   |
| chr3 (178936091) | PIK3CA | p.E545K              | 9.0%        | 7.00 – 10.60%    |
| chr3 (178952085) | PIK3CA | p.H1047R             | 17.5%       | 14.00 – 21.00%   |
| chr10 (43613843) | RET    | p.L769L              | 60.0%       | 54.00 – 66.00%   |
| chr17 (7579472)  | TP53   | p.P72R               | 92.5%       | 83.25 – 100.00%  |

## Material Provided

Mimix™ OncoSpan FFPE Reference Standard

- Catalogue/Model number: HD832-IVD  
One vial, 1x FFPE curl, 15 µm thickness.



## Precautions and Warning

- This product is not classified as dangerous, however; users are advised to handle all materials as potentially biohazardous, following standard laboratory safety procedures.
- Dispose of all waste materials in accordance with local regulations.
- Avoid contamination of the product when opening and closing the vial by making sure that the curl is intact and sits at the bottom of the vial.
- Inspect the vial for any signs of melting before opening.

## Storage and Handling

- Store unopened at 2 to 8 °C.
- Do not use this product beyond the expiration date printed on the product label.
- Please see the following instructions for handling:
  - Open the tube by twisting and lifting the lid.
  - If needed, please remove the curl carefully without destruction by using suitable forceps (e.g. anatomical forceps) and transfer the curl appropriately.
  - If the whole curl is to be removed from the vial, we recommend all handling takes place at temperatures below room temperature (~15-25 °C).
  - If the curl does not need to be transferred, please keep it in the vial and perform your assay accordingly (e.g. gDNA extraction).

**Quality Control and Performance Characteristics**

Variant detection when using the Mimix™ OncoSpan FFPE Standard may vary depending on the library preparation method, sequencing platform, and bioinformatics pipeline used for NGS assays, as well as primer/probe set, analysis type/platform, and assay conditions for ddPCR. The expected variant list of the OncoSpan FFPE provided can be used to compare each subsequent run.

Following our QC procedure, the acceptance criteria are described below (Table 2). Users should note that results will vary depending on the equipment and workflow used. Users are expected to perform their own validation to determine AFs within the workflow to be monitored. The below reported data were obtained using the validated equipment and workflow as specified for each quality control test.

Table 2. List of quality control tests applied and the corresponding acceptance criteria.

| Test name             | Test purpose                             | Acceptance criteria                                 |
|-----------------------|--|---|
| Fluorometric analysis | Measure quantity and quality of FFPE DNA | dsDNA ≥400ng  |
| ddPCR                 | Measure stability of AF                  | Within the accepted target range. Refer to Table 1. |

The gDNA yield obtained from the OncoSpan FFPE Reference Standard is shown in Figure 1.

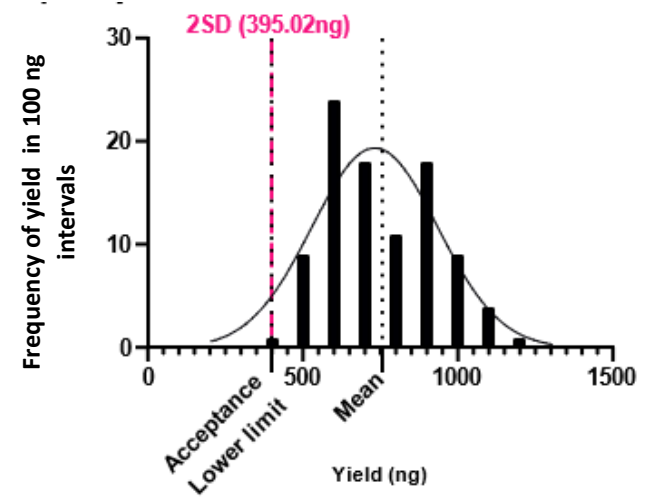


Figure 1. Frequency distribution for gDNA amount extracted from OncoSpan FFPE Reference Standard. The mean and lower acceptance criteria are shown.

Batch to batch variability of gDNA yield extracted from OncoSpan FFPE Reference Standard is shown in Figure 2.

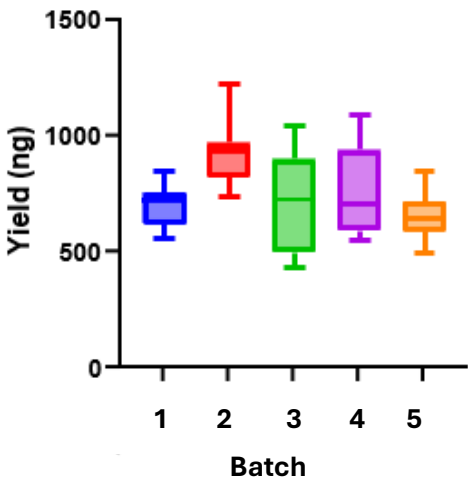


Figure 2. Batch-to-batch yield of gDNA extracted from OncoSpan FFPE Reference Standard.

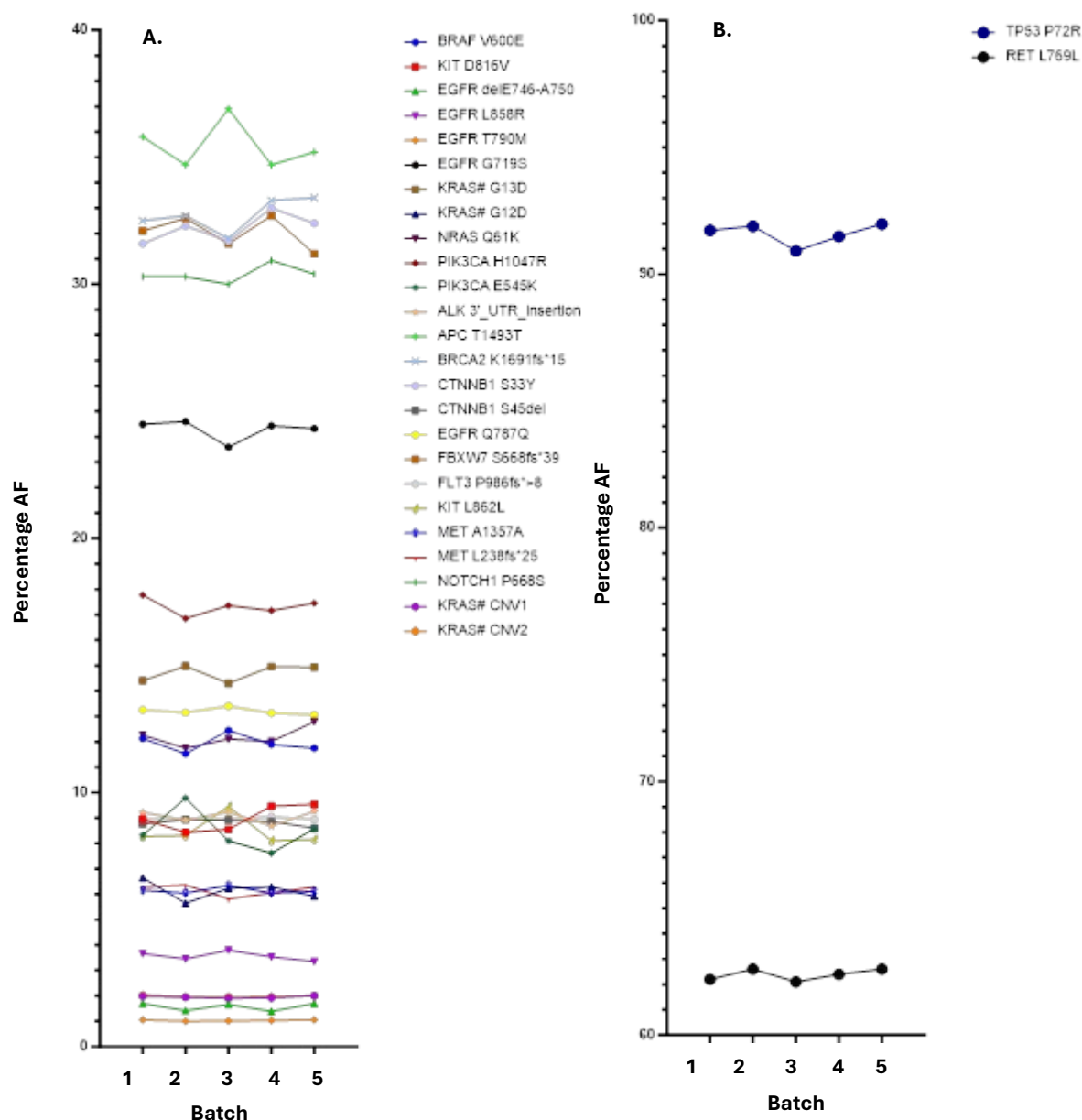


Figure 3. (A and B) Batch to batch variation of AF for different genes and mutations in the OncoSpan gDNA Reference Standard confirmed by ddPCR.

#### Limitations

- The Mimix™ OncoSpan FFPE Reference Standard is not to be used as a substitute for the internal controls provided by manufacturers within their IVD assay kits.

#### Technical Support

Scientific support

Tel: 800-235-9880 (Option 2) Or 303-604-9499 (Option 2)

Fax: 1-800-292-6088 Or 303-604-9680

Email: [technical@horizondiscovery.com](mailto:technical@horizondiscovery.com)

#### Customer support

Tel: 800-235-9880 (Option 1) Or 303-604-9499 (Option 1)

Fax: 1-800-292-6088 Or 303-604-9680

Email: [orders@horizondiscovery.com](mailto:orders@horizondiscovery.com)

For a high-confidence variant list go to <https://horizondiscovery.com/reference-standards/products/oncospan-ffpe-reference-standard-ivd>

Batch-specific whole exome sequencing data is provided upon purchase as an online link.

References

1 Vinkšiel, M., Writzl, K., Maver, A. *et al.* Improving diagnostics of rare genetic diseases with NGS approaches. *J Community Genet* 12, 247–256 (2021). <https://doi.org/10.1007/s12687-020-00500-5>


2 Boycott K, Hartley T, Adam S, et al. The clinical application of genome-wide sequencing for monogenic diseases in Canada: Position Statement of the Canadian College of Medical Geneticists. *J Med Genet.* 2015;52(7):431-437. doi:10.1136/jmedgenet-2015-103144


3 Mirabile A, Sangiorgio G, Bonacci PG, et al. Advancing Pathogen Identification: The Role of Digital PCR in Enhancing Diagnostic Power in Different Settings. *Diagnostics (Basel).* 2024;14(15):1598. Published 2024 Jul 25. doi:10.3390/diagnostics14151598


Glossary


- IVD


In Vitro Diagnostic Medical Device
- REF

Catalogue Number
- 

Manufacturer
- 

Temperature Limit
- 

Caution
- 

Consult instructions for use
- 

Use-by date
- UDI

Unique device identifier
- QTY

Quantity
- CONTROL+

Positive control

Limited are indirectly wholly owned subsidiaries of Revvity, Inc.

Mimix and OncoSpan are trademarks of Revvity Health Sciences Inc.



Revvity Discovery Limited  
8100 Cambridge  
Research Park, Waterbeach,  
Cambridge, CB25 9TL  
United Kingdom

Document ID: 03-09-204  
Version: 00

**Trademark and Patent information**

©2025 Revvity Health Sciences Inc. All rights reserved.  
Horizon Discovery Biosciences Limited is a wholly owned subsidiary of Revvity Discovery Limited. Revvity Discovery Limited and Horizon Discovery Biosciences