

# Spotlight Myeloid™ Liquid Biopsy Panel

## *Comprehensive and exon-level hotspot coverage of 23 clinically relevant myeloid genes*

The Spotlight Myeloid Panel combines content from peer-reviewed publications and other sources to offer comprehensive and exon-level hotspot coverage of 23 clinically-relevant genes implicated in acute myeloid leukemia (AML), myelodysplastic syndrome (MDS) and myeloproliferative neoplasms (MPN). Because of the high prevalence of myeloid variants in clonal hematopoiesis of indeterminate potential (CHIP) studies, this is an ideal panel for uncovering CH variants in either cell-derived or cell-free DNA samples.

**Comprehensive:** Covers over 12,000 COSMIC and 1,900 ClinVar mutations and generates targeted libraries compatible with Illumina® sequencing platforms.

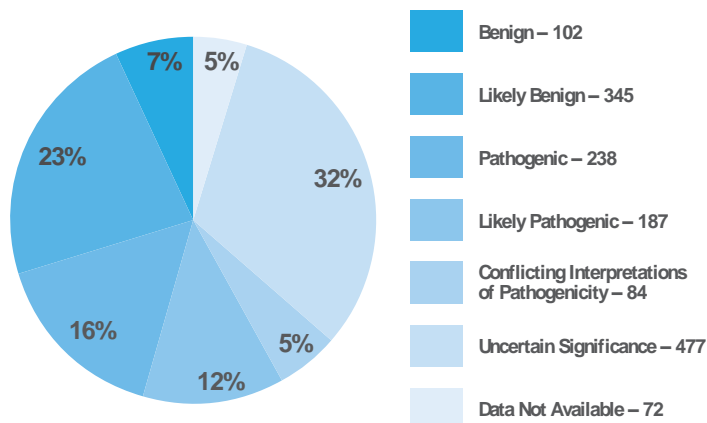
**Informative:** Provides full exon coverage of TP53 with flanking intron/exon boundaries.

**Powerful:** Simultaneous detection of single nucleotide variants (SNV) and small insertion and deletions (indels), all within a fast and easy single-tube assay workflow.

**Flexible:** Add your own targets using our pre-validated primers or others content. The possibilities are endless.

**Other key features include:**

- Robust performance on cell-free DNA and FFPE samples starting with as little as 20 ng input.
- Highly sensitive detection of rare variants down to 0.1% allele frequency.
- Sequence ready libraries in just 2 hours.
- Streamlined, free analysis via Fluxion’s ERASE-Seq Variant Caller.
- Compatible with all Illumina platforms.



**Figure 1.** ACMG and AMP-classified variants of clinical significance covered by the Myeloid Panel.

**This panel enables:**

- Full exon coverage of TP53
- Compatibility with a wide range of samples
- Free data analysis using the proprietary ERASE-Seq caller
- Addition of your own custom content

## Product Specifications

| Feature                           | Specification   |
|-----------------------------------|---|
| Input DNA Required                | 20 ng   |
| Starting Sample                   | cfDNA, buffy coat, FFPE   |
| Library Prep Time Required        | 2 hours   |
| Number of Amplicons               | 478   |
| Amplicon Size                     | Average 142 bp  |
| Total Target Size                 | 42 kb   |
| Number of Genes                   | 23  |
| <b>Gene List</b>                  | ASXL1, CALR, CEBPA, CSF3R, DNMT3A, EZH2, FLT3, HRAS, IDH1, IDH2, RUNX1, SETBP1, SF3B1, SRSF2, JAK2, JAK3, KDM6A, KIT, MPL, NPM1, RUNX1, SETBP1, SF3B1, SRSF2, TET2, <b>TP53</b> , U2AF1 |
|                                   | (Bold indicates whole CDS coverage)   |
| Sensitivity                       | 0.1% AF and below   |
| Validated Sequencers              | MiSeq, HiSeq, NextSeq, Iseq, NovaSeq  |
| Average Sequencing Depth Required | 15,000X per sample  |

## Ordering Information

| Description   | Part Number |
|---|-------------|
| 24-sample Spotlight Myeloid Panel, including primer pool, library kit, and adapter kit; 0.1% AF sensitivity, 20 ng input, 2 reaction tubes per sample, cloud-based ERASE-Seq variant calling included | 910-0133    |

For research use only. Not for use in diagnostic procedures.