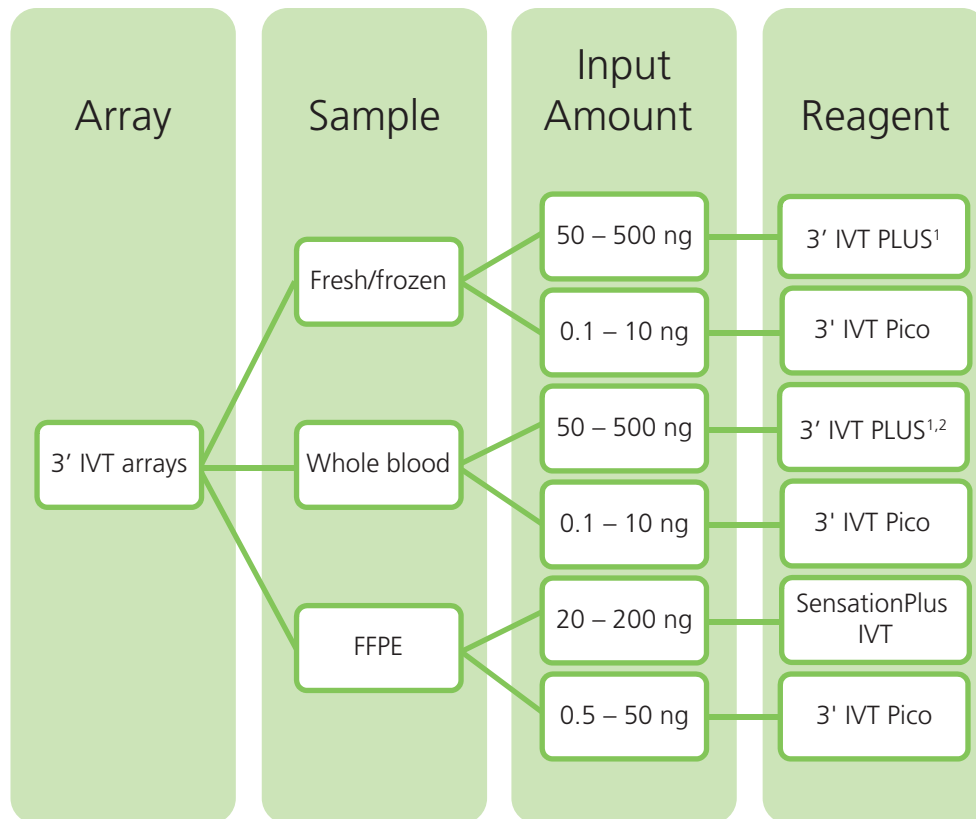


Expression microarray reagent guide

3' IVT assays



GeneChip® 3' IVT PLUS Kit (3' IVT PLUS)¹

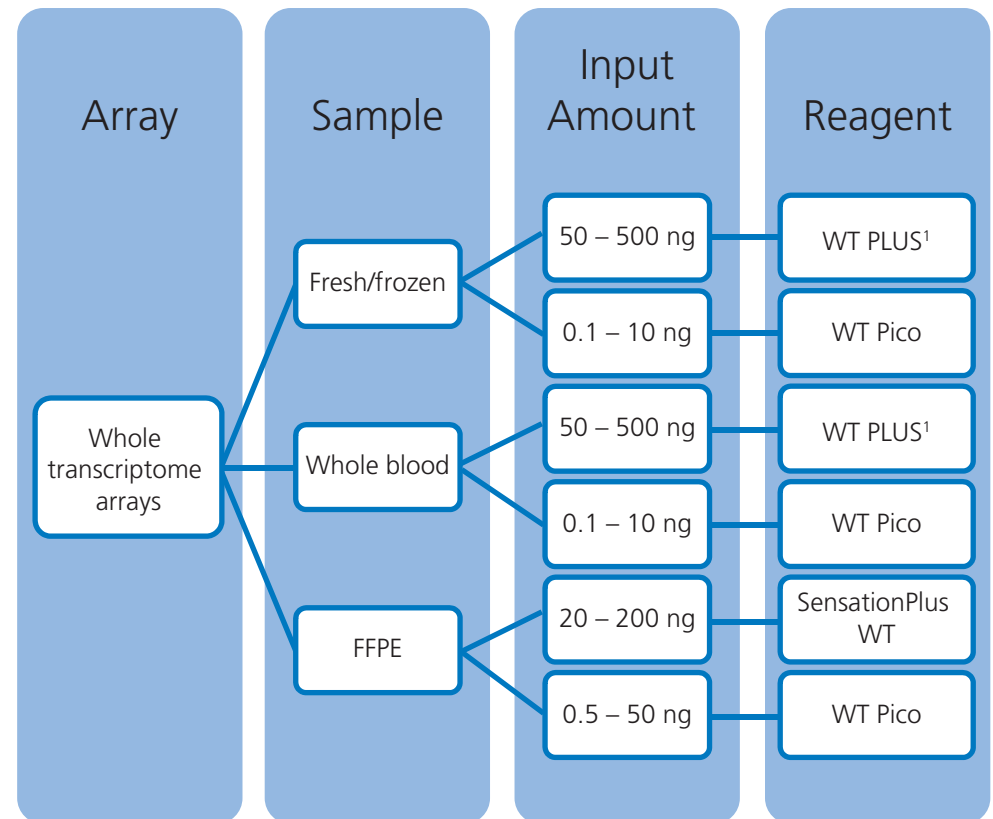
SensationPlus™ FFPE Amplification and 3' IVT Labeling Kit (SensationPlus IVT)

GeneChip® 3' IVT Pico Kit (3' IVT Pico)

¹Also available as a high-throughput (HT) format for automation.

²Requires globin mRNA reduction.

Whole transcriptome assays



GeneChip® WT PLUS Kit (WT PLUS)¹

GeneChip® WT Pico Kit (WT Pico)

SensationPlus™ FFPE Amplification and WT Labeling Kit (SensationPlus WT)

For Research Use Only. Not for use in diagnostic procedures.



Expression microarray reagent guide

miRNA assays

Array	Sample	Input Amount	Reagent
miRNA arrays	Fresh/frozen	100 – 1000 ng	FlashTag
	Whole Blood	100 – 1000 ng	FlashTag
	FFPE	100 – 1000 ng	FlashTag

Affymetrix FlashTag™ Biotin HSR RNA Labeling Kits (FlashTag)

Microarray categories*

Whole-transcriptome (WT) arrays

GeneChip® Human Transcriptome Array 2.0 (HTA 2.0)
 GeneChip® mouse and rat transcriptome arrays 1.0 (MTA 1.0, RTA 1.0)
 GeneChip® human, mouse, and rat gene 2.0 ST arrays
 GeneChip® human, mouse, and rat gene 1.0 ST arrays
 Model and applied research organisms gene 1.0 ST arrays
 GeneChip® human, mouse, and rat exon 1.0 ST arrays

3' IVT arrays

GeneChip® Human Genome U133 Plus 2.0 Array
 GeneChip® PrimeView® Human Gene Expression Array
 GeneChip® Mouse Genome 430 2.0 Array
 GeneChip® Rat Genome 230 2.0 Array
 Model and applied research organism genome arrays

miRNA arrays

GeneChip® miRNA 4.0 Array
 GeneChip® miRNA 3.0 Array
 GeneChip® miRNA 2.0 Array

*List includes most common array types only and is not comprehensive. Find additional WT and 3' IVT arrays as well as information about the MyGeneChip™ Custom Array Program at www.affymetrix.com.