

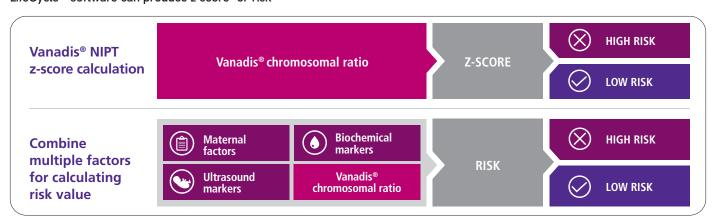
Flexible. Reliable. LifeCycle.

LifeCycle[™] prenatal screening software

Ideal solution for all prenatal screening laboratories.

Flexible screening	 Multiple screening strategies are supported Software is able to combine the demographics, biomarker data and Vanadis chromosomal ratios to produce risks. 			
Tailored to population	Optimize risk calculation to your populationAbility to level or change MoM equations			
Accessible	Accessible via web browser or over remote connection outside laboratory			
Configurable	High level of configurability and connectivity options			

LifeCycle™ software can produce z-score* or risk

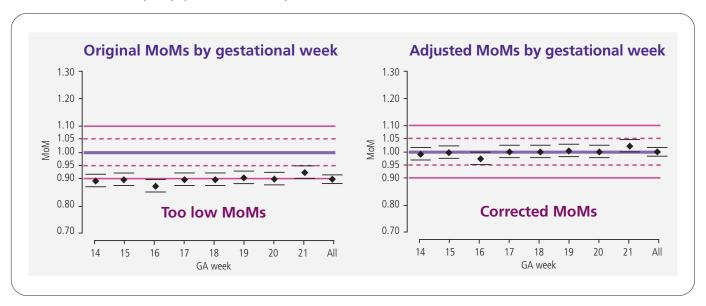


^{*}Vanadis® NIPT support from version 7.0 onwards

Products may not be licensed in all countries such as the United States.

Please check with your local representative for availability.

Tailor risk calculation to your population with LifeCycle™ software



LifeCycle™ software versions

Feature	v. 6.0	v. 7.0	v. 7.0 rev 2	v. 7.0 rev 3
User-configurable	✓	✓	✓	✓
Browser-based	✓	✓	✓	✓
Remote users	✓	✓	✓	✓
User-configurable file based LIMS import/export	✓	✓	✓	✓
T21, T18, T13 & pre-eclampsia risk	✓	✓	✓	✓
What if feature	✓	✓	✓	✓
Optional HL7 import / export for LIMS and Ultrasound equipment	✓	✓	✓	✓
Z-score calculation for Vanadis® NIPT system		✓	✓	✓
Fetal sex determination for Vanadis® NIPT system*		✓	✓	✓
Retest screening strategy for Vanadis® NIPT system			✓	✓
Ability to print labels for blood tubes			✓	✓
Pre-eclampsia twin risk calculation			✓	✓
sFlt-1/PlGF ratio calculation				✓

^{*}Sex determination is offered as an option for nonmedical purposes, only in countries where permitted. Sex determination is not offered in India under PNDT regulation.



