

RELEASE NOTES

November 15, 2017

Precision ID NGS System

Precision ID GlobalFiler NGS STR Panel v2 and Analysis Files
Converge™ Software 2.0 – Kinship and NGS Data Analysis Modules

TABLE OF CONTENTS

- Products Affected
- Software Overview
- New Features
- Installation/Upgrade
- Known Issues and Limitations
- System Requirements
- Release & Compatibility Summary
- Legal Notices and Disclaimer

PRODUCTS AFFECTED

- A33114 Precision ID GlobalFiler NGS STR Panel v2 Analysis Files
- Converge Software 2.0
 - Kinship and Paternity Analysis Module
 - Next-Generation Sequencing Data Analysis Module

SOFTWARE OVERVIEW

Converge Software is a modular, enterprise platform that provides key analysis to perform next-generation sequencing (NGS) and complex kinship and paternity testing. The software serves as a central repository for case data and simplifies the daily complexities of forensic case management.

Converge NGS Data Analysis module is required to generate profiles from the Applied Biosystems Precision ID GlobalFiler NGS STR Panel v2. NGS analysis functionality includes information on Short Tandem Repeats (STRs) including allele calls, sequence motifs, Single Nucleotide Polymorphisms (SNPs) in flanking regions. NGS data uses Applied Biosystems GeneMapper ID-X Software inspired process quality values (PQVs) and flags such as allele number (AN), off-ladder allele (OL), peak height ratio (PHR), below stochastic threshold (BST), and control concordance (CC). Preconfigured analysis settings are provided and may be modified by the laboratory as needed. Additionally, full auditing functionality is included for chain of custody requirements. Using the NGS Data Analysis module with the Case Management application allows for simplified comparison of NGS and CE profiles.

For Research Forensic or Paternity Use Only. Not for use in diagnostic procedures. For licensing and limited use restrictions visit thermofisher.com/HIDlicensing.

Converge Kinship and Paternity module integrates with Converge Software and GeneMapper ID-X genotyping software to help enable automated analysis, reporting of routine paternity and other relationship testing cases, and allows for genetic likelihood ratio calculation, as well as reporting.

Case Management application supports case, subject, genotype profile, and laboratory data management, and allows for automated data transfer and integration with various forensic DNA laboratory systems.

NEW FEATURES

Modifications have been made to Precision ID GlobalFiler™ NGS STR Panel v2 and its analysis files.

INSTALLATION/UPGRADE

TSS

It is recommended that your Torrent Server, Ion Chef, and Ion S5/S5 XL remain on TSS 5.2.2. For TSS v5.2.2 support documentation, refer to the NGS Software Support page [here](#).

BED/JSON Files

The following files must be installed onto TSS v5.2.2 before analyzing data generated with Precision ID GlobalFiler NGS STR Panel v2:

The following files have been updated to v1.1 (previously v1.0):

- Precision_ID_GlobalFiler_NGS_STR_Panel_Target_v1.1.bed
- Precision_ID_GlobalFiler_NGS_STR_Panel_Hotspot_v1.1.bed
- Precision_ID_GlobalFiler_NGS_STR_Panel_AnalysisParams_v1.1.json
- Precision_ID_GlobalFiler_NGS_STR_Control_Sample_male007_v1.1.json (Template-1002 in Converge)
- Precision_ID_GlobalFiler_NGS_STR_Control_Sample_9947A_v1.1.json (Template-1003 in Converge)
- Precision_ID_GlobalFiler_NGS_STR_Control_Sample_NegCtrl_v1.1 (Template-1004 in Converge)
- Precision_ID_GlobalFiler_NGS_STR_Control_Sample_9947A_and_male007_and_NegCtrl_v1.1 (Template-1005 in Converge)

Download the analysis files [here](#).

Once the zip file has been downloaded, it can be unzipped and installed on the Torrent Suite server.

For Research Forensic or Paternity Use Only. Not for use in diagnostic procedures. For licensing and limited use restrictions visit [thermofisher.com/HIDlicensing](https://www.thermofisher.com/HIDlicensing).

Under the “Admin” tab in Converge, in the “Add-ons” section, in the “Manage Add-ons” sub-section, please disregard the following files and refer to the newest versions provided in links.

- Precision_ID_GlobalFiler_NGS_STR_Panel_Target_v1.0
- Precision_ID_GlobalFiler_NGS_STR_Panel_Hotspot_v1.0
- Precision_ID_GlobalFiler_NGS_STR_Panel_AnalysisParams_v1.0

Under the “Admin” tab in Converge, in the “Global Settings” section, in the “Manage Templates” sub- section, please disregard the following templates and use the new versions of the files instead.

- Template-1002, Template-1003, Template-1004, Template-1005

KNOWN ISSUES AND LIMITATIONS

Converge Software 2.0

1. Microsoft™ Excel files exported from the STR grid in the results page do not filter alleles showing zero coverage. The results page in the user interface accurately filters these instances.

23	D5S818	10, 13	PASS	PASS	PASS	PASS	Sequence	Long Sequence
	Allele		Status	Coverage				
	10		ABOVE_ST	3008			[AGAT]10	D5S818[CE10]-chr5-hg19 123111250-123111293 [AGAT]10
	10			8			[AGAT]10	D5S818[CE10]-chr5-hg19 123111250-123111293 [AGAT]10
	10			3990			[AGAT]10	D5S818[CE10]-chr5-hg19 123111250-123111293 [AGAT]10 123111306-G
	10			2			[AGAT]10	D5S818[CE10]-chr5-hg19 123111250-123111293 [AGAT]10 123111306-
	10			5			[AGAT]10	D5S818[CE10]-chr5-hg19 123111250-123111293 [AGAT]10 123111306-G 123111309-
	10			2			[AGAT]10	D5S818[CE10]-chr5-hg19 123111250-123111293 [AGAT]10 123111214- 123111306-G
	13		ABOVE_ST	2057			[AGAT]13	D5S818[CE13]-chr5-hg19 123111250-123111293 [AGAT]13
	13			6			[AGAT]13	D5S818[CE13]-chr5-hg19 123111250-123111293 [AGAT]13
	13			2046			[AGAT]13	D5S818[CE13]-chr5-hg19 123111250-123111293 [AGAT]13 123111306-G
	13			2			[AGAT]13	D5S818[CE13]-chr5-hg19 123111250-123111293 [AGAT]13 123111305- 123111306-G
	13			3			[AGAT]13	D5S818[CE13]-chr5-hg19 123111250-123111293 [AGAT]13 123111306-
	13			6			[AGAT]13	D5S818[CE13]-chr5-hg19 123111250-123111293 [AGAT]13 123111306-G 123111309-

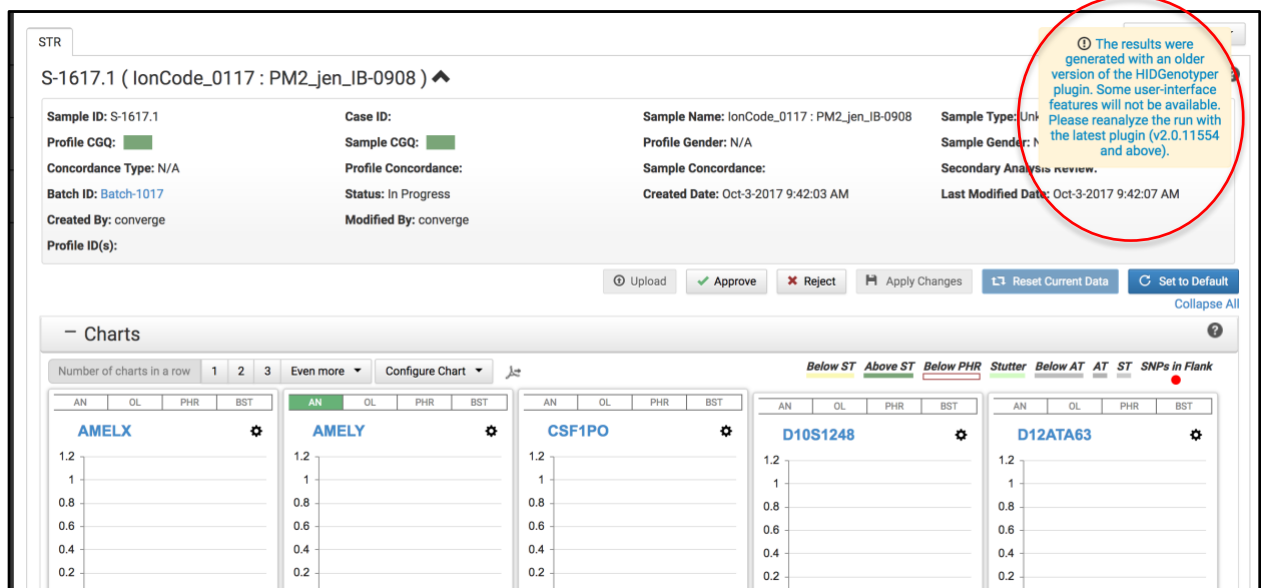
2. Excel files exported from the STR grid in the results page display some alleles with coverage percentage as ‘N/A’. This occurs when the forward strand coverage of alleles is zero. The results page in the user interface has the correct coverage percentage displayed.

2.3.0	Long Sequence	Ref/Alt	Read	Coverage	% Quality
2.3.1	Penta_D[CE10]-chr21-hg19 45056086-45056150 [TCCTT]10			N/A	
2.3.2	Penta_D[CE13]-chr21-hg19 45056086-45056150 [TCCTT]13			N/A	

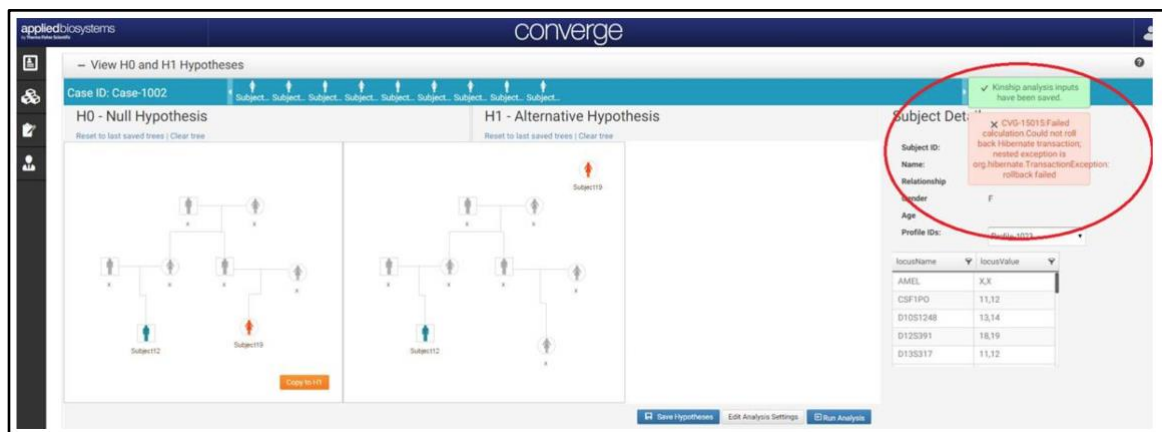
3. When a sample has no data after secondary analysis, Converge user interface displays the following message:

‘The results were generated with an older version of the HIDGenotyper plugin. Some user interface features will not be available. Please analyze the run with the latest plugin (v2.0.11554 and above).’

This error message is incorrect and should be ignored.



- Transaction time out error might occur when using the kinship module for analysis that may exceed 1 minute.



For Research Forensic or Paternity Use Only. Not for use in diagnostic procedures. For licensing and limited use restrictions visit thermofisher.com/HIDlicensing.

- CGQ flag does not account for the result of the concordance check. When the concordance (profile, sample level) has failed, indicated by red flag, on the results page on Converge, the CGQ (profile/sample level) is green, it should be red instead.

STR Select Sample

NTC-2017 (IonCode_0102 : NTC)

Sample ID: NTC-2017	Case ID:	Sample Name: IonCode_0102 : NTC	Sample Type: Negative Control
Profile CGQ: ■	Sample CGQ: ■	Profile Gender: No Male Presence	Sample Gender: No Male Presence
Concordance Type: NTC	Profile Concordance: ■	Sample Concordance: ■	Secondary Analysis Review:
Batch ID: Batch-1073	Status: In Progress	Created Date: Sep-21-2017 10:38:07 AM	Last Modified Date: Sep-21-2017 10:38:14 AM
Created By: converge	Modified By: converge		
Profile ID(s):			

- Under the “Admin” tab in Converge, in the “Add-ons” section, in the “Manage Add-ons” sub-section, please disregard the following: target, hotspot and analysis parameters files. To obtain these, refer to the URL provided in the installation/upgrade section above.

Admin

Kinship and Paternity

- Loci Inclusion List
- Analysis Parameters Settings
- Manage Population Database

Global Settings

- Manage Licenses
- Manage Loci
- Manage Kits
- Manage Templates
- Manage Organizations
- Manage Custom Fields
- Manage Case Archives
- Configure Search

Security

- Manage User Accounts
- Manage User Roles
- Manage User Account Security Policies

Add-ons

- Manage Add-ons

Manage Add-ons

More

Page 1 of 1

10

Target System	Name	Description	Size
TSS v5.2.2	HIDGenotyper.zip	TSS Plugin v2.0	333.8 MB
TSS v5.2.2	Precision_ID_GlobalFiler_N...	PrecisionID GlobalFiler NGS STR Panel Hotspot	4.3 KB
TSS v5.2.2	Precision_ID_GlobalFiler_N...	PrecisionID GlobalFiler NGS STR Panel Targets	3.6 KB
TSS v5.2.2	Precision_ID_GlobalFiler_N...	PrecisionID GlobalFiler NGS STR Analysis Parameters	14.8 KB
TSS v5.2.2	Control_Samples.zip	Control Samples 2.0	2.5 KB

For Research Forensic or Paternity Use Only. Not for use in diagnostic procedures. For licensing and limited use restrictions visit thermofisher.com/HIDlicensing.

SYSTEM REQUIREMENTS

TSS

- TSS v5.2.2

Converge Software Server

- Dell™ PowerEdge™
- Red Hat™ Enterprise Linux™ operating system
- Apache™ Tomcat™ application server that runs on Converge software v2.0
- PostgreSQL database server that stores the data for the server and software
- Google™ Chrome™ browser
- Automatic configuration of IP, domain name service (DNS), and Windows internet name service (WINS) settings via dynamic host configuration protocol (DHCP)

Supported Browsers

- Microsoft™ Internet Explorer™ v11
- Google Chrome™ 55
- Apple™ Safari™ v8

Recommended Software (not provided)

- Adobe™ Acrobat Reader
- Microsoft Excel

RELEASE & COMPATIBILITY SUMMARY

Updates in **BOLD** below.

SYSTEM	TYPE	DESCRIPTION	CURRENT VERSION	PREVIOUS VERSION
TSS	Software	TSS Compatibility	v5.2.2	N/A
Genotyper	Plugin	HIDGenotyper	v2.0_r16121	N/A
Control Publisher	Publisher	Control_Samples	v1.01	N/A
	Control File	Precision_ID_GlobalFiler_NGS_ST R_Control_Sample_male007	v1.1	N/A
	Control File	Precision_ID_GlobalFiler_NGS_ST R_Control_Sample_9947A	v1.1	N/A
	Control File	Precision_ID_GlobalFiler_NGS_ST R_Control_Sample_NegCtrl	v1.1	N/A
	Control File	Precision_ID_GlobalFiler_NGS_ST R_Control_Sample_9947A_and_male007_and_NegCtrl	v1.1	N/A
BED/ JSON File	Target	Precision_ID_GlobalFiler_NGS_ST R_Panel_Target.bed	v1.1	v1.0
	Hotspot	Precision_ID_GlobalFiler_NGS_ST R_Panel_Hotspot.bed	v1.1	v1.0
	Analysis Parameters	Precision_ID_GlobalFiler_NGS_ST R_Panel_AnalysisParams.json	v1.1	v1.0
Converge	Platform	Platform	Serial No. 16151	N/A
	License	Platform	2.0	1.0
	Module	Case Management	2.0	2.0
	Module	Kinship and Paternity Analysis	1.0	1.0
	License	Kinship	1.0	1.0
	Module	Next-Generation Sequencing	1.0	N/A
	License	NGS	1.0	N/A

For Research Forensic or Paternity Use Only. Not for use in diagnostic procedures. For licensing and limited use restrictions visit thermofisher.com/HIDlicensing.

LEGAL NOTICES AND DISCLAIMER

For Research, Forensic, or Paternity Use Only. Not for use in diagnostic procedures. For licensing and limited use restrictions visit thermofisher.com/HIDlicensing.

Information in this document is subject to change without notice.

DISCLAIMER: LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE (S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NONINFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

TRADEMARKS

© 2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.