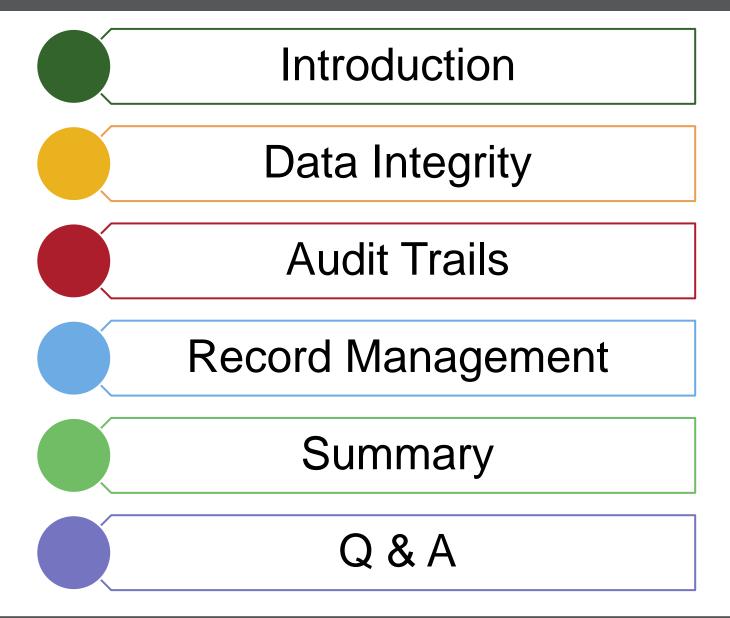


ThermoFisher SCIENTIFIC

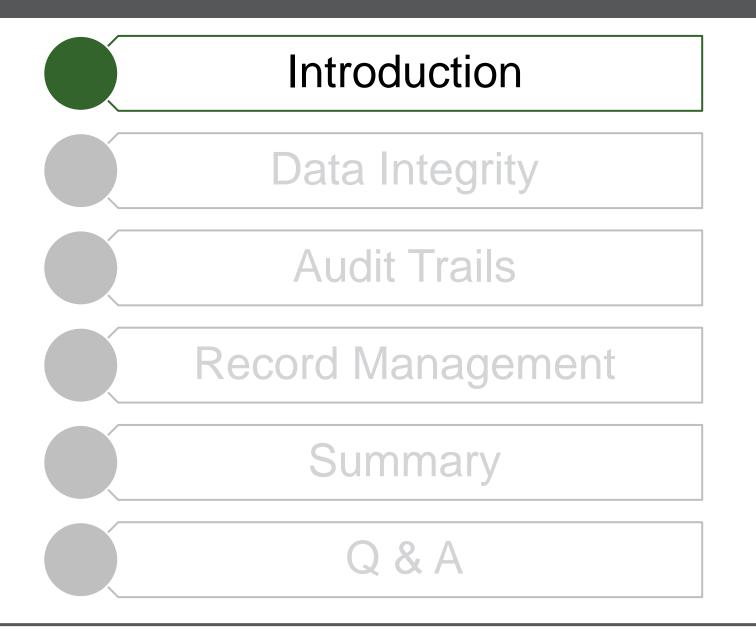
Comply with Data Integrity Regulations with Chromeleon CDS Software

Darren Barrington-Light Product Marketing Specialist Chromatography Software

Content Overview



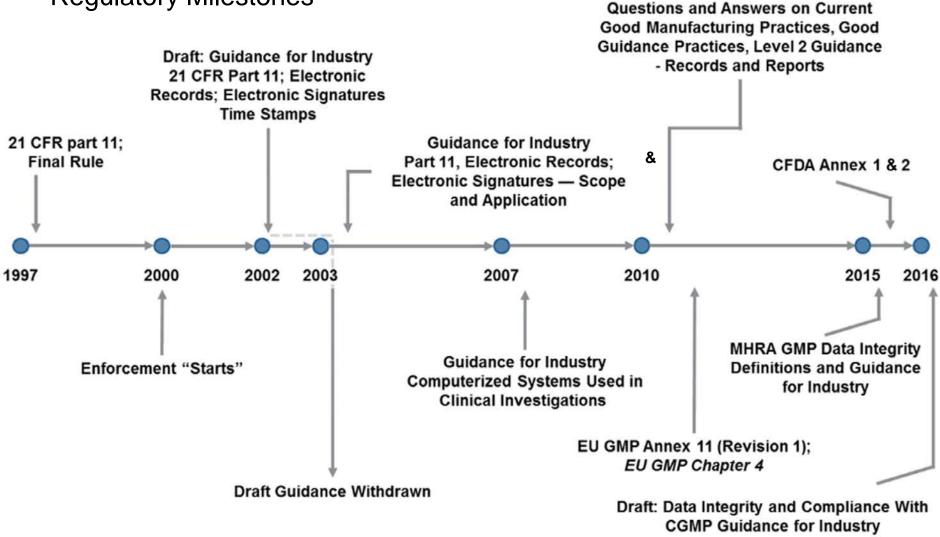






Introduction

Regulatory Milestones



Data integrity requirements

.S. Food and Drug Administration

- New guidelines in 2015 & 2016 from MHRA, CFDA and FDA
- Covers authenticity and security of data
- Clarifies role of data integrity in current good manufacturing practice (CGMP) for drugs
- Contains current thinking on creation and handling of data in accordance with CGMP requirements
- Software needs to meet regulatory requirements.
 - Includes requirements for 21 CFR Part 11 compliance

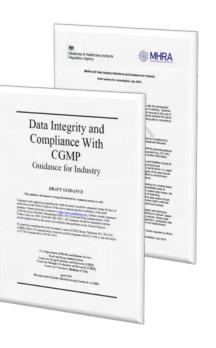








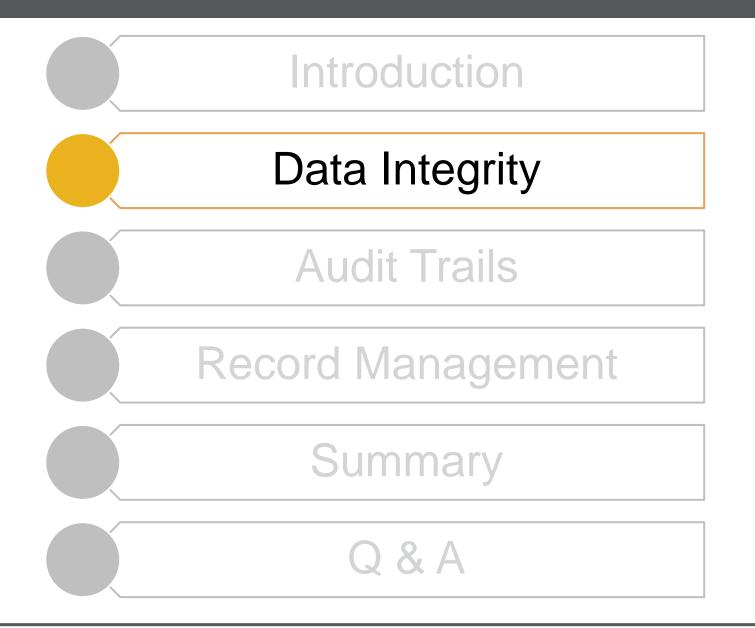






- In recent years, FDA and MHRA have increased focus on data integrity during CGMP inspections.
- Observed CGMP violations involving data integrity during inspections
- Ensuring data integrity is important
 - Industry's responsibility to ensure safety, efficacy, and quality of drugs
 - Regulatory responsibility to protect the public health
- How can Thermo Scientific[™] Chromeleon[™] Chromatography Data System (CDS) software help?







Data Integrity

- What is "data integrity"?
 - Completeness, consistency, and accuracy of data

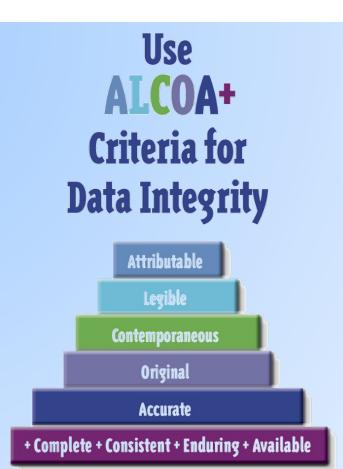
Attributable Who performed an action & when?

Legible Can you read the data file & written entries throughout the life cycle?

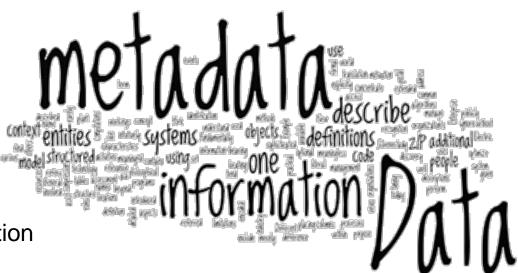
Contemporaneous Documented at the time of the activity.

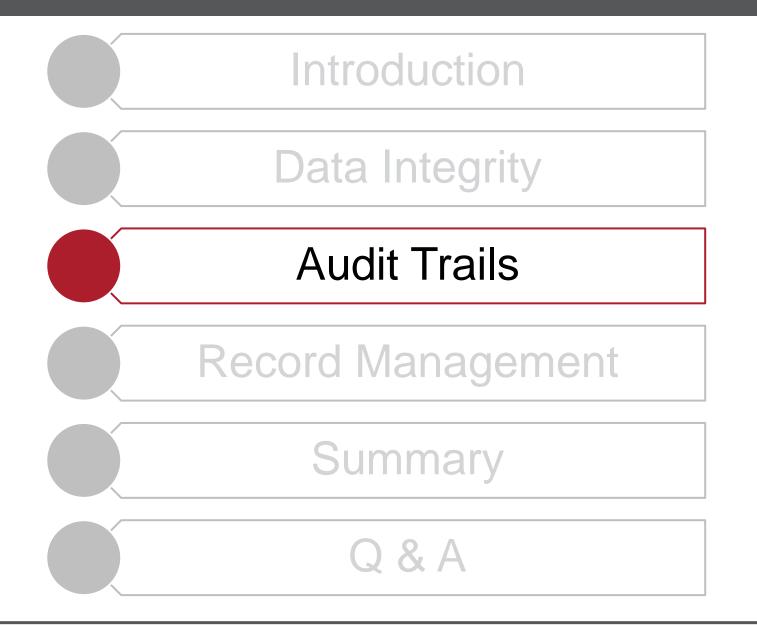
Original Original record or a certified copy

Accurate No errors or editing without documented amendments



- What is "data"?
 - Any recorded information. 2 types:
 - Static fixed-data document such as paper record or electronic image
 - **Dynamic** record format that allows interaction between user and record content
 - E.g. chromatogram allowing user to change baseline or reprocess data so peaks may change size; user modifiable formulas or entries in spreadsheet
- What is "metadata"?
 - Metadata is data about data
 - Contextual information required to understand data
 - A data value is meaningless by itself without additional information
 - Metadata can describe a single piece of data, a data set or collection

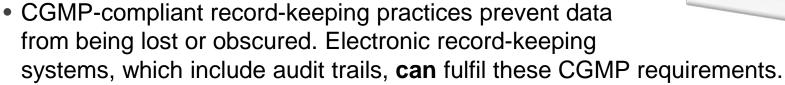






Audit Trails

- What is an "audit trail"?
 - Secure, computer-generated, time-stamped electronic record that allows for reconstruction of events relating to creation, modification, or deletion of an electronic record.
 - Electronic audit trails include those that track creation, modification, or deletion of data **and** those that track actions at the record or system level.



Chronology of the "who, what, when, and why" of a record



- Is it acceptable to only save final results from reprocessed laboratory chromatography?
 - No. Analytical methods should be capable and stable
 - For most lab analyses, reprocessing data should not be regularly needed
 - If chromatography is reprocessed, written procedures must be established and followed and each result retained for review
 - Must have complete data in lab records, which includes raw data, graphs, charts, and spectra from lab instruments
- Predicated rules show this refers to reprocessing of **samples**, not **data**.
 - Must keep all data (and metadata) relating to each analysis
 - Versioning helps track changes to objects



Chromeleon CDS provides industry-leading audit trails

- Detailed audit trails for:
 - Chromeleon Console & Chromatography Studio
 - Administration Console, Instrument Configuration Manager, Data Vault Manager
 - Chromeleon Station
- Logs <u>all</u> user, instrument & system actions
 - E.g. user logon/logoff, instrument setup, sequence start/stop, data processing, admin changes, services startup/shutdown, etc.

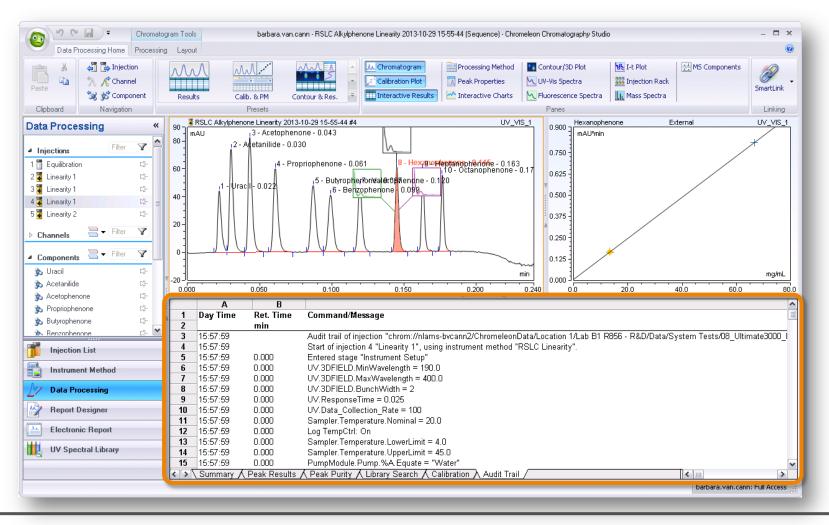
	Report 😥 Show Ch	nanges 🦃 R	estore l	nclude Sul	ofolde	rs from 1	2/09/2016	🔲 🔻 to 10/	11/2016 🖉 🍸	Filtering 📇 Grou	ping	+		
ĺ	Name	Туре	Version	Da	ite / Tir	ne 👻	Operator	Operation	Comment	Additio	nal Information			
Î	🚰 ChromeleonLocal	Data Vault	1	10/11/20	16 12:3	2:43 +00:00	cmadmin	Changed	Lockout other users	Access control cl	hanged			
	ChromeleonLocal	Data Vault	1	10/11/20	16 10:0	9:57 +00:00	cmadmin	Changed		Enable versioning	g and data audit tr	ails,		
	Instrument Utilization	Folder	1	12/09/20	16 14:0	4:39 +01:00	cmadmin	Created		Imported from C:	Users\darrenbl\D	eskto		
						Cate	Jory	Da	ate/Time 🔺	Operator	Host Name	Object	Operation	
					U	ser Manag	ement	10/11/201	6 09:53:46 +00:00	cmadmin		Full Access	🌚 Changed Role	
	_	-			▶ Lo	gon and S	ignature	10/11/201	6 09:54:04 +00:00	cmadmin		cmadmin	ogoff	
					Lo	gon and S	ignature	10/11/201	6 09:54:14 +00:00	cmadmin		cmadmin	🕜 Logon	
								. <u>.</u>		<u>i</u>	<u>i</u>	£		



- Chromeleon CDS consolidated audit trail viewer
 - Combines all Audit Trails within Organizational Unit
 - Events are categorized for filtering and grouping
 - Entries contain significant detail

🕉 Chromeleon Administration Co	onsole (UKALT-DBL-SRFP4)							
License Manager (localhost)	🗄 🔁 🔜 <u>R</u> eport 💯 Sh	ow Changes <u>F</u> rom 09/11/2	016 🔲 🔻	<u>T</u> o <u>11/11/201</u>	16 🔲 🔻	Y Filtering 🔤 Groupir	ng	
Organizational Units	Category	Date/Time 🔺	Operator	Host Name	Object	Operation	Comment	Additi
⊡	Logon and Signature	09/11/2016 16:43:12 +00:00	cmadmin		cmadmin	of Logon		Logon Role =
⊕ 🥵 User Database () eWorkflow Tags	Logon and Signature	09/11/2016 16:45:19 +00:00	cmadmin		cmadmin	- J Logon		
Audit Trail	Global Policies	09/11/2016 16:45:48 +00:00	cmadmin		<global></global>	Changed		Changed Poli
⊞⊶∰ QA ⊞⊶∰ R&D	User Management	09/11/2016 16:46:36 +00:00	cmadmin		<global></global>	Changed User		Changed Poli
Domain Resources	Logon and Signature	09/11/2016 16:48:04 +00:00	cmadmin		cmadmin	Logon		Logon Role =
Computers	Logon and Signature	09/11/2016 16:49:54 +00:00	cmadmin		cmadmin	Logoff		
- Instruments	Logon and Signature	09/11/2016 16:50:02 +00:00	cmadmin		cmadmin	Jogon		Logon Role =
Chromeleon Consoles	Logon and Signature	09/11/2016 17:07:24 +00:00	cmadmin		cmadmin	Logoff		
	Logon and Signature	10/11/2016 08:33:58 +00:00	cmadmin		cmadmin	👸 Logon Failed		Count of wron
	Logon and Signature	10/11/2016 09:31:30 +00:00	cmadmin		cmadmin	✓ Logon		Count of wron

• Instrument daily event audit trail for each injection



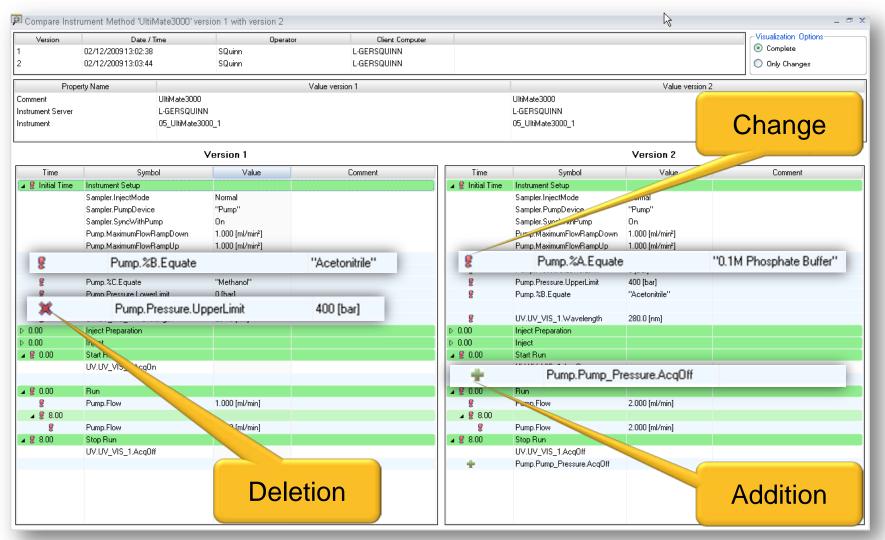


- Modification history audit trail
- Displays complete change history of each object (versions)
- Can roll back (restore) a previous version

a 1	ta Audit Trail - Diuretics Method Scouti	0		er 🏆 Clear Filters 🗮 Grouping	(#A	a, a		0
±	Name	Туре	Version	Date / Time	Operator	Operation	Comment	1
17 🗂	Sequence Diuretics Method Scouting	Sequence	169	25/11/2010 01:31:01 PM +01:00	CmUserAms	Changed Injection (3)		1
17 📝	SCOUTING	Processing Method	168	25/11/2010 11:46:12 AM +01:00	CmUserAms	Copied	From 'chrom://cm_bvc/Chromeleo	
17 /		Processing Method	167	25/11/2010 11:46:12 AM +01:00	CmUserAms	Deleted		1
17 🎁	Sequence Diuretics Method Scouting	Sequence	166	23/11/2010 05:08:51 PM +01:00	CmUserAms	Copied	From 'chrom://c-ams-cm7/Central	٦
17 🍸	Sequence Diuretics Method Scouting	Sequence	166	23/11/2010 04:57:16 PM +01:00	CmUserAms	Deleted Signal (302)		1
17 🎁	Sequence Diuretics Method Scouting	Sequence	166	23/11/2010 04:47:45 PM +01:00	CmUserAms	Exported	Sent to: C:\Documents and Settin	
17 🍸	Sequence Diuretics Method Scouting	Sequence	166	23/11/2010 04:07:42 PM +01:00	CmUserAms	Changed	Changed	1
17 🎁	👖 Sequence Diuretics Method Scouting	Sequence	165	23/11/2010 03:52:31 PM +01:00	CmUserAms	Changed	Changed	1
17 /	SCOUTING	Processing Method	164	23/11/2010 03:35:26 PM +01:00	CmUserAms	Changed		
16 🎁	Sequence Diuretics Method Scouting	Sequence	163	23/11/2010 03:10:00 PM +01:00	CmUserAms	Renamed Injection (29)	From 'Diuretics' to 'Method 1'.	
16 🎁	Sequence Diuretics Method Scouting	Sequence	162	23/11/2010 02:15:51 PM +01:00	CmUserAms	Changed Injection (29)		
16 🙀	🌠 Default	View Settings	161	23/11/2010 01:59:57 PM +01:00	CmUserAms	Changed		
16 🎁	Sequence Diuretics Method Scouting	Sequence	160	23/11/2010 01:49:51 PM +01:00	CmUserAms	Changed		
16 🔛	Default DAD	Report Template	160	23/11/2010 01:49:51 PM +01:00	CmUserAms	Created		
16 /	7	Processing Method	159	23/11/2010 01:07:07 PM +01:00	CmUserAms	Changed		
16 🔨	⊵ 8390.smp/UV_VIS_3.channel/UV_VIS_3	Chromatogram	158	23/11/2010 01:04:03 PM +01:00	CmUserAms	Changed		
16 🛝	ע 8390.smp/UV_VIS_3.channel/UV_VIS_3	Chromatogram	157	23/11/2010 12:18:56 PM +01:00	CmUserAms	Changed		
16 🔨	20653.smp/UV_VIS_3.channel/UV_VIS_3	Chromatogram	156	23/11/2010 12:18:16 PM +01:00	CmUserAms	Changed		
-	⊵ 8482.smp/UV_VIS_3.channel/UV_VIS_3	Chromatogram	155	23/11/2010 12:17:40 PM +01:00	CmUserAms	Changed		
15 🛆	└ 11926.smp/UV_VIS_3.channel/UV_VIS_3	Chromatogram	154	23/11/2010 12:17:25 PM +01:00	CmUserAms	Changed		
15 🛆	26757.smp/UV_VIS_3.channel/UV_VIS_3	Chromatogram	153	23/11/2010 12:16:52 PM +01:00	CmUserAms	Changed		
15 🛆	ע 5998.smp/UV_VIS_3.channel/UV_VIS_3	Chromatogram	152	23/11/2010 12:16:24 PM +01:00	CmUserAms	Changed		
15 🛆	22383.smp/UV_VIS_3.channel/UV_VIS_3	Chromatogram	151	23/11/2010 12:15:13 PM +01:00	CmUserAms	Changed		
	21095.smp/UV_VIS_3.channel/UV_VIS_3	Chromatogram	150	23/11/2010 12:14:52 PM +01:00	CmUserAms	Changed		
-	28723.smp/UV_VIS_3.channel/UV_VIS_3	Chromatogram	149	23/11/2010 12:13:08 PM +01:00	CmUserAms	Changed		
15 🛆	└ 1634.smp/UV_VIS_3.channel/UV_VIS_3	Chromatogram	148	23/11/2010 12:12:18 PM +01:00	CmUserAms	Changed		
15 🔥	28723.smp/UV_VIS_3.channel/UV_VIS_3	Chromatogram	147	23/11/2010 12:11:46 PM +01:00	CmUserAms	Changed		



Version comparison tool





• File section in injection list and folder view shows current version

🔮 cmadmin - Chromeleon Console				- 🗆	×					
😋 Back 💿 <u>C</u> reate <u>F</u> ile <u>E</u> dit <u>V</u> iew <u>T</u> ools <u>H</u> elp					0					
Data «	Runs									
Image: Station Qual 2016-11- Image: Station Qual 2016-11-	Type Date Modified 10 Sequence 10/11/2016 09:49:51 +00:00	Version	Commer tation Operation	nt onal Qualification						
È- È- StationQualification È- È- Runs ▼	🔮 cmadmin - Chromeleon Console								- 🗆	×
Station Qual 2016-11-10 User training	Sack 🔘 Create File Edit V	iew <u>T</u> ools <u>H</u> elp								(
P Deleted Items	Data «				Station Qual 20	16-11-10				
	Filter 🝸	Run Finished					<u></u>	CM_OQ>	(Not Connecte	ed)
Instruments	ChromeleonLocal	🗄 🚽 Save 🎯 Studio 🖨) Print 👻 🖻	Up 🛛 📴 Insert	t Row 👻 🏨 Fill Down	Lock	🝸 Filtering 📒	Grouping		;
Data	Analytics Report Templates Difference Instrument Data	# UV_VIS_1 ▶ N	lame	1	Туре	Level	Position	Volume [µl]	Instrument Metho	od H
eWorkflows	. My Sequences		Alkylphen	ones 1	Calibration Standard	01	RA1	2.000	CM_OQ	Rack View
	🖻 🧰 Runs	2	Alkylphen	ones 2	Calibration Standard	01	RA1	4.000	CM_OQ	2
Folder 'Runs' selected	User training	3	Alkylphen	ones 3	Calibration Standard	01	RA1	8.000	CM_OQ	
	Deleted Items	4 A	Alkylphen	ones 4	Calibration Standard	01	RA1	12.000	CM_OQ	
		5	Alkylphen	ones 5	Calibration Standard	01	RA1	16.000	CM_OQ	
		6 M M	Alkylphen	ones 6	Calibration Standard	01	RA1	20.000	CM_OQ	4
		7 ΛΛΛ	Alkylphen	ones 7	Unknown		RA2	4.000	CM_OQ	
					Click here to add a n	ew injection		•	.•	
		•)	▶ «
	Instruments	Name		Туре	Date Me	⊲ odified	Version	С	omment	-
	Data			Processing Met Instrument Meth	·····		1			
	eWorkflows			Report Template View Settings	e 10/11/2016 09:49 10/11/2016 09:49		1			
		Associated Items	Custom Se	quence Variables						
	Sequence 'Station Qual 2016-11-10' selecte	ed							:madmin: Full Ac	ccess

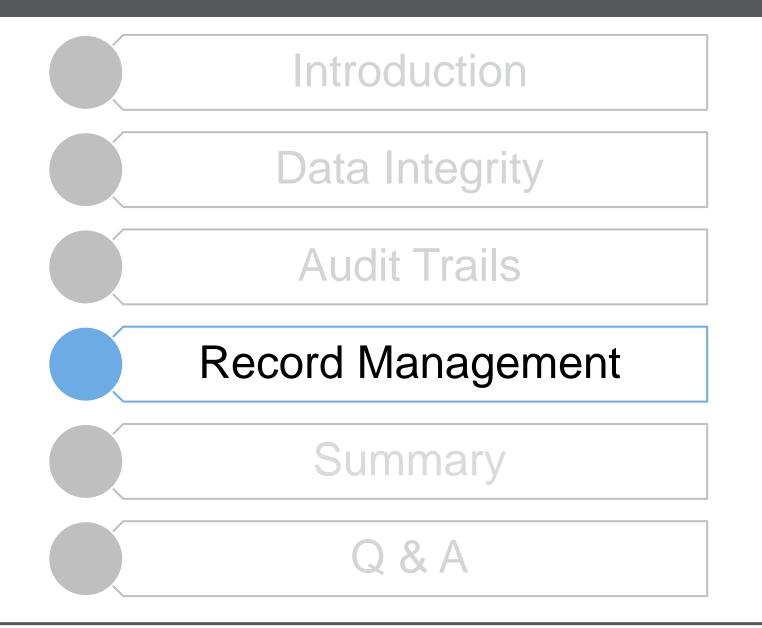


Privileged Actions

- Require authorization for specified actions
- Ensure users add appropriate comments for specified actions
- Use standardized comments or free entry

🕉 Chromeleon Administration Co	onsole (UKALT-DBL-SRFP4)		- 0	×
Gobal Policies Gobal Policies Gobal Chromeleon Users	Save Policies Report Report		al Units	0
Access Groups	Lipection Locking Remote Data Vaults	Settings Standard Comments	J Authorization and Comment Required	×
Power User TI Admin To Admin	 Instrument Audit Trails Detection Algorithm Version User Templates Location Watermark UI Customizations Privileged Actions Email Configuration Station Audit Raw Data 	Action Instrument Configuration Instrument Method Create Delete Modify Copy	With the second seco	
		Move Rename Instrument Audit Trail Data Vault Folder		221 characters remaining







All data should be stored, controlled and archived correctly

- When does electronic data become a CGMP record?
 - When generated to satisfy any CGMP requirement, all data becomes a CGMP record
 - Must document, or save, data at time of performance (contemporaneous)
 - E.g. chromatograms should be sent to long-term storage upon run completion
 - Not acceptable to record data on paper that will be discarded after data are transcribed to permanent laboratory notebook
 - Not acceptable to store data electronically in temporary memory, in manner that allows for manipulation, before creating permanent record

All data should be stored, controlled and archived correctly

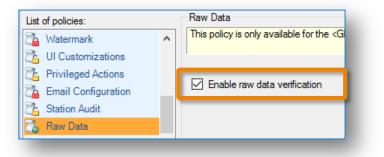
- How do regulatory bodies use the term "backup"?
 - A true copy of original data that is maintained securely throughout records retention period.
 - Backup file should contain data (with associated metadata) in original format or in format compatible with original format.
- Must not be confused with backup copies!
 - Created during normal computer use and temporarily maintained for disaster recovery (e.g. in case of a computer crash or other interruption)
- How does Chromeleon CDS help?

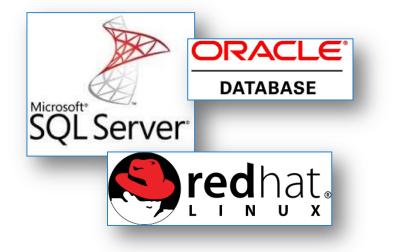




Immediate transfer to long-term storage

- Chromeleon CDS uses relational database
 - Microsoft[®] SQL[®] Server Express (2005 SP3, 2008 R2 SP2 & 2014 SP2)
 - Microsoft[®] SQL[®] Server (2008 R2, 2012 & 2014 SP2)
 - Oracle[®] Database 11g & 12c
 - Oracle[®] Database 12c on Red Hat[®] Enterprise Linux[®] 7
- Detect changes to raw files
 - Automatic detection of raw files being manipulated outside Chromeleon CDS

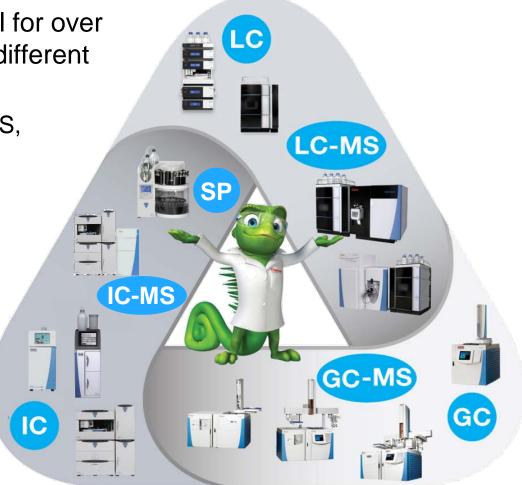




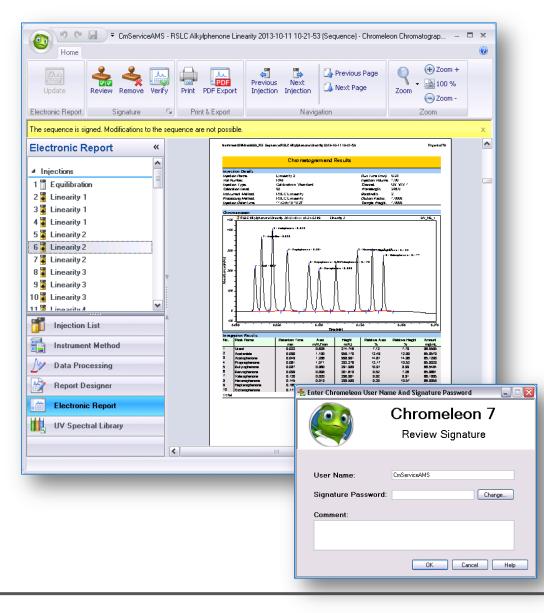


Record Management – Data Storage

- Bi-directional instrument control for over 450 different modules from 17 different manufacturers including:
 - Thermo Scientific (LC, IC, GC, MS, Sample Prep)
 - Agilent (LC, GC)
 - Waters (LC)
 - Shimadzu (LC, GC)
 - Varian (LC, GC)
 - Perkin Elmer (LC, GC)
 - Gilson (LC)
 - CTC PAL (Sample Prep)
- Highest data transport security
 - Communication secured using latest standards



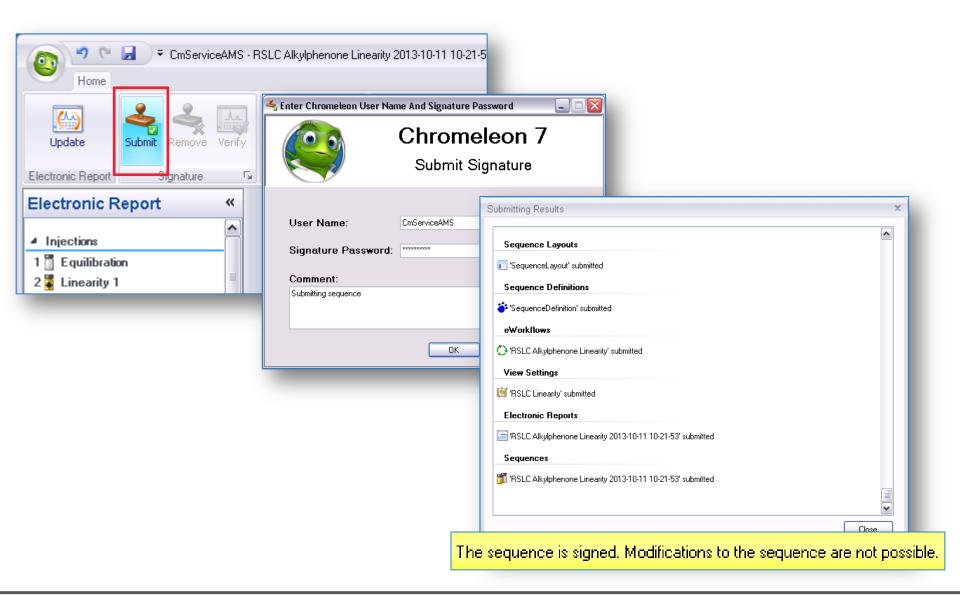
Record Management – Data Control



- Electronic Report
 - Snapshot in time of sequence
 - Any changes tracked
 - Can be sent for review
- Electronic Signatures
 - 3 Levels of signature Submit, Review, Approve
 - Applied to electronic report
 - Complete sequence protected by checksums
 - Controlled by privileges



Record Management – Data Control

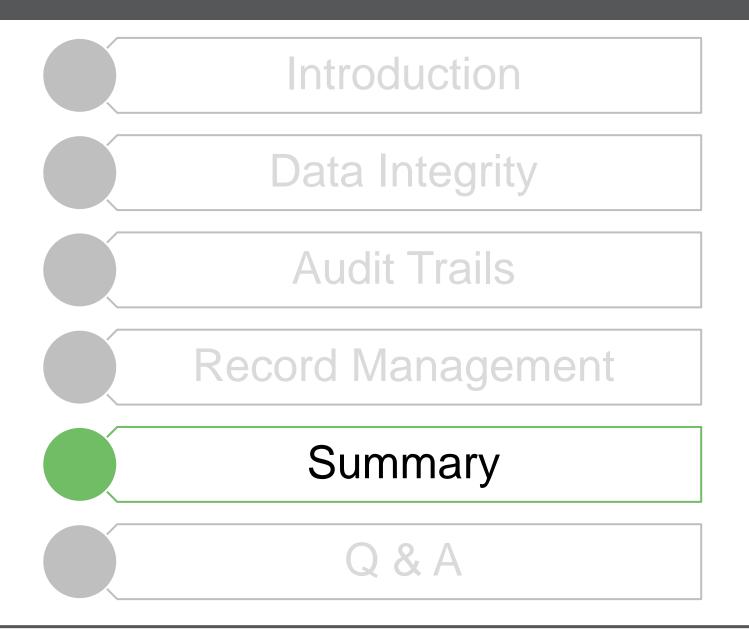




- Chromeleon Scheduler automatically archive/backup data using powerful Injection Queries
 - Transfer to another Data Vault direct access with full metadata
- Export to cmbx or other format (e.g. PDF, GAML, XLS, etc.)

Schromeleon Administration Console (n	nlams-cm02)	- = X	
E-G License Manager	Scheduler is running.	Last retrieval of service data: 23-Dec-13 14:43:52	
	Task List Audit Trail Service Account Service Log Abo	out	
Scheduler Scheduler (Chromeleon Domain)	2 Tasks Enabled Task Name Status	Chromeleon Administration Console (nl	 ams-cm02)
Garage Concerns and the concerns of the c	Archive Data older than one month Waiting	🖃 🗔 License Manager	Task is done. Next start: 23-Jan-14 14:26:24 Last retrieval of task data: 23-Dec-13 44:44:57 @
	Cleanup cached network sequences Waiting	Client Licenses	General Schedule Items Action Audit Trail
I Full Access			Name:
Operator Cm Administrator		Archive Data older than one mor	Archive Data older than one month
	1	User Database Orromeleon Users	Owner: Barbara van Cann
Testing			Description: Archive all data that is older than one month to the central archiving data vault.
Global Policies		Full Access	
Local Machine		Operator Cm Administrator	
		Laborant	C Enabled
		🔤 😳 Testing 🔤 🚛 Audit Trail	Specify additional settings that affect the behavior of the task.
		Global Policies	✓ If the task fails, restart every: 1 hour
			Attempt restart up to: 3 🛟 times
			Terminate Stop Start Delete Discard changes Save changes



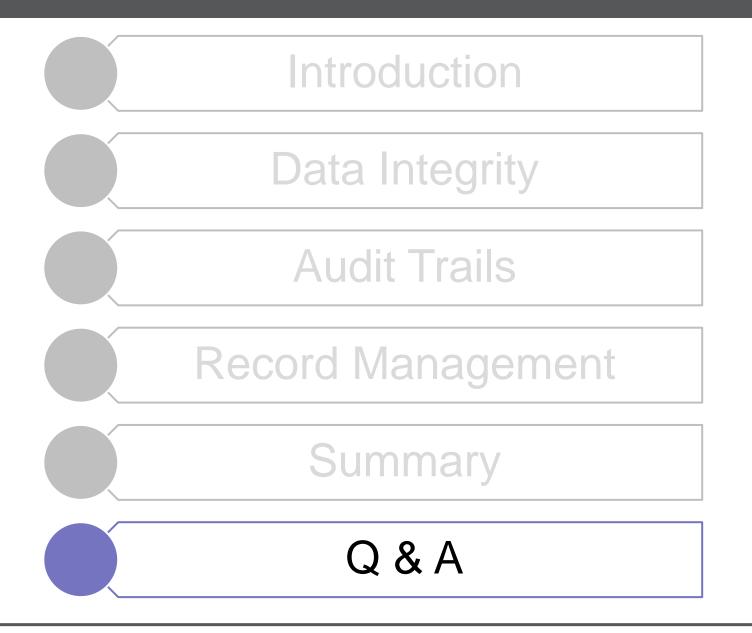




Summary

- Data integrity and compliance go hand-in-hand
 - Subject to elevated scrutiny by regulatory bodies
- Chromeleon CDS meets all of your data integrity needs:
 - Ensures completeness, consistency, and accuracy of data
 - ALCOA+
 - Comprehensive audit trails
 - Who, what, when and why of data
 - Maintains record authenticity
 - Contemporaneous collection
 - Automated long-term data retention with easy accessibility







For more information about Chromeleon 7.2 CDS or to request an informational 'demo' CD, visit: *thermofisher.com/Chromeleon*

Like Charlie Chromeleon on Facebook to follow him and get important updates on chromatography software!



facebook.com/CharlieLovesChromatography



Any questions?





Do you have additional questions or do you want to talk to an expert from Thermo Fisher Scientific?

Please send an E-Mail to analyze.eu@thermofisher.com and we will get back to you.

