

NIR Technology at Freeport McMoRan Mine Sites

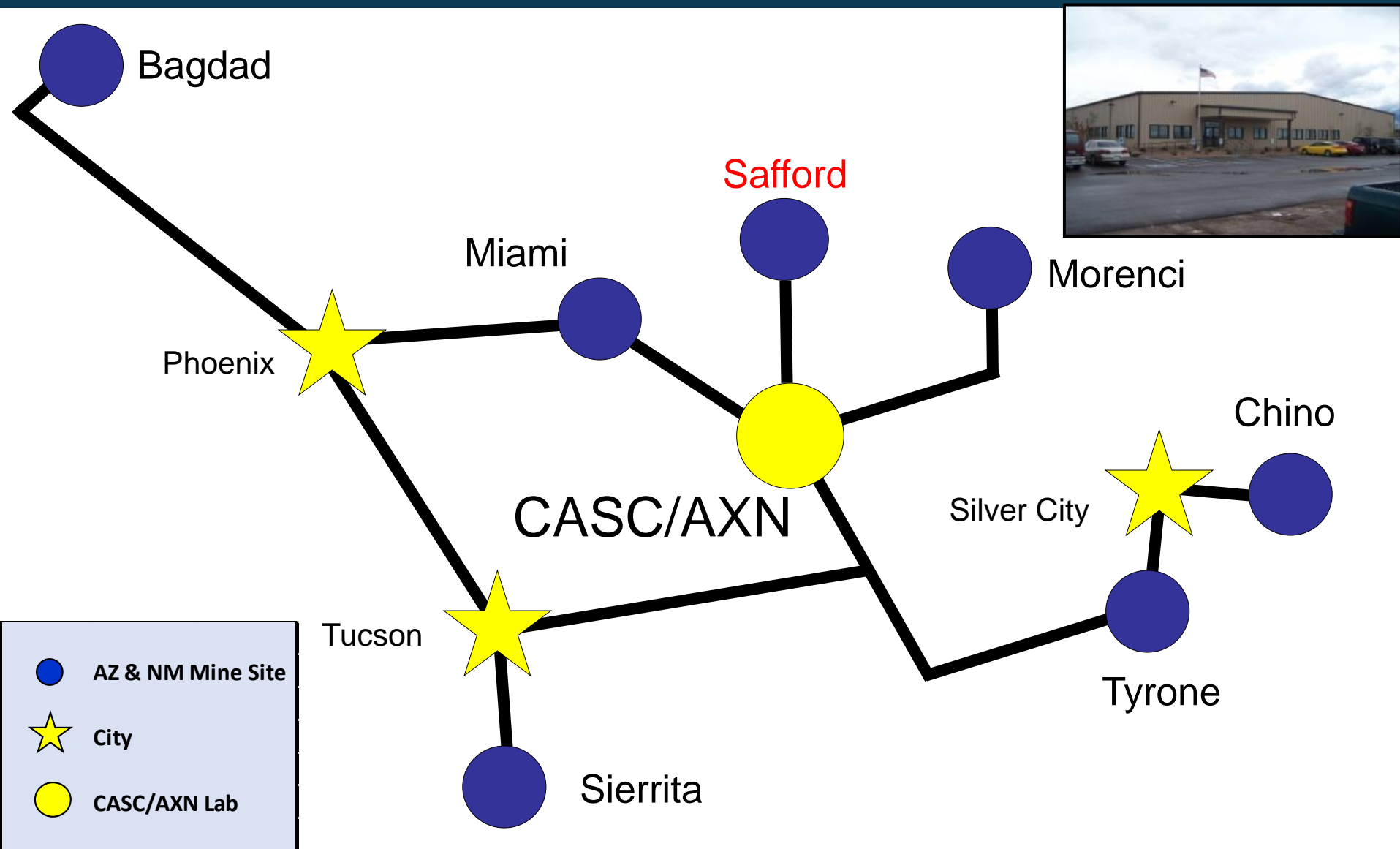
***Julie Schaub
July 21, 2020***

- Overview of Freeport McMoRan's Mineralogy Department
 - Central Analytical Service Center (CASC)
 - Automated X-ray & NIR (AXN)
 - Technology Center Tucson (TCT)
- Technology Comparisons
- QA/QC
- NIR Data Use at Morenci and Safford mine sites

Overview

North American Operations Served by CASC and AXN

FREEPORT
FOREMOST IN COPPER



Overview of Freeport Technology Center Mineralogy Department

AXN

- X-ray Diffraction Analysis (XRD)
- Near Infrared Analysis (NIR)
- Cation Exchange Capacity (CEC)

Process Mineralogy Lab

- Optical Microscopy
- Scanning Electron Microscopy
- Automated Mineralogy



Automated X-ray / NIR Overview (AXN)

- Primary Instrumentation:
 - 7 X-ray diffraction (XRD) analyzers
 - 4 Near Infrared (NIR) analyzers
 - 3 Ion Chromatographs (CEC: Cation Exchange Capacity)
 - Automated pulverization circuit
- Provides bulk, clay and alteration mineralogy
- High throughput/fast turnaround/highly automated
- Linked to CASC robotic lines for daily blast-hole analyses
- Turn Around Time as fast as <24 hours
- Average workload of 3,000 XRD and 16,000 NIR analyses per month



AXN Robotic Circuit

Technology Comparisons

Factors

- Cost
- TAT
- Difficulty of Analysis

XRD/NIR - rapid, cost-effective non-destructive, accurate, & efficient analytical method making it ideal for linking with CASC robotic lines for daily blasthole analysis

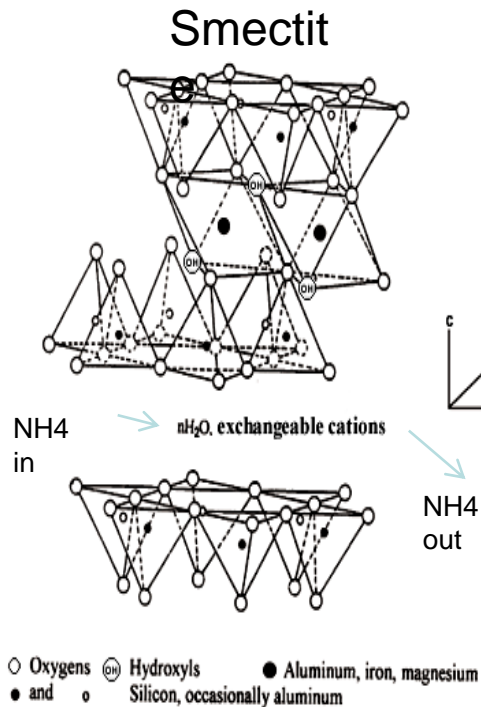
QEMSCAN provides valuable information not obtained through XRD/NIR, but requires more complex sample preparation and analysis

- Labor intensive sample preparation, slower TAT
- Costly
- Detection limits under normal operating parameters
 - Unable to detect approx. <8 Wt.% elemental Cu in minerals
 - Requires SEM analysis to better define SIP definitions
 - Light element limitations (oxygen, carbon, hydrogen)
- Beam resolution on fine textures & phase intergrowths

- Detection limits vary by mineral. Typically $\leq 0.5\%$
 - Low grade copper minerals fall below detection limits
 - Not an ideal tool for Cu deportment
- Minor phases overlapped by major peaks - resolved by reconciliation with assay
- Preferred orientation of micas/molybdenite during press mounting
 - Press mounts are back filled
 - Topas software allows calculation for correction

- Unable to detect amorphous non-crystalline material
 - Can be sample prep generated or natural
 - Use of CEC to estimate swelling capacity of sample (swelling clay)
 - Micronization - control amorphous & reduce microabsorption effects

- Not applicable for Cu mineral department
- Limited by reference data – XRD & CEC
- Requires site specific calibration model & periodic model updates for quantitative predictions
- Limited spectrum, limited constituent set (clays, micas, chlorite)
- Reference data required for calibration models : XRD and CEC



- Measures the quantity of cations (Mg, K, Ca, & Na) that a sample can exchange.
- Predicts swelling capacity and reported as swelling clay content.
- Most clays tough to quantify via XRD (amorphous).
- Normalized CEC value to XRD data for bulk mineralogy.
- Final measurements by ion chromatography.



Standard Mineral CEC Values

Kaolinite	3-15	
Halloysite 2H ₂ O		5-10
Halloysite 4H ₂ O		40-50
Montmorillonite-group	70-100	
Illite	10-40	
Vermiculite	100-150	
Chlorite	10-40	
Glauconite	11-20+	
Palygorskite-group	20-30	
Allophane	~70	

QAQC

- NIST Corundum standard analyzed daily on each XRD
 - EVA peak position and area plotted on control chart
- AXN Ore Standard (Granite) analyzed daily on each XRD
 - Synthetic standard being developed for XRD
 - Topas Rietveld refinement results for charting
- Minimum of 5% of samples are duplicated
- Blank (quartz sand) samples prepared daily to check for contamination
- Pulverizing mills checked weekly to ensure proper particle size is produced
- New development - Reconciliation of Rietveld analysis with chemical assay

- Mylar Standard analyzed daily on NIR
- Site Specific Ore Standards analyzed daily on NIR
- M-distance >5 = Failed Model built into LIMS
 - Failed Model samples get CEC
- Random samples selected for XRD/CEC; Quarterly XRD/CEC vs. NIR correlation reports issued
- Model updates performed based on XRD-CEC/NIR correlations & excessive Failed Models

Data Use

- NIR data used daily for shovel routing and ore blending & routing (critical)
- XRD data
 - Retroactive requests in areas that are identified “problematic”. Helps the mill to prepare ahead of time when they get back to these areas
 - Requests randomly selected to understand full bulk mineralogy in all areas

- Daily blast hole submittals for NIR and select XRD
 - NIR – Basic gangue for rock type
 - CEC/NIR Clay – blending
 - XRD – Bulk and alteration mineralogy
- NIR, XRD, and CEC (clay) data used in geologic model
 - Model is provided to the engineers for ore routing, blending, and forecasting
 - Data is referenced at least once a day
- Minerals identified include: swelling clay, muscovite, biotite, kaolinite, chlorite
- NIR is quick & inexpensive (\$6/sample), but requires spectral model to be built based upon XRD & CEC data
- Requires site specific calibration model & periodic model updates for quantitative predictions

Ore Control Modeling

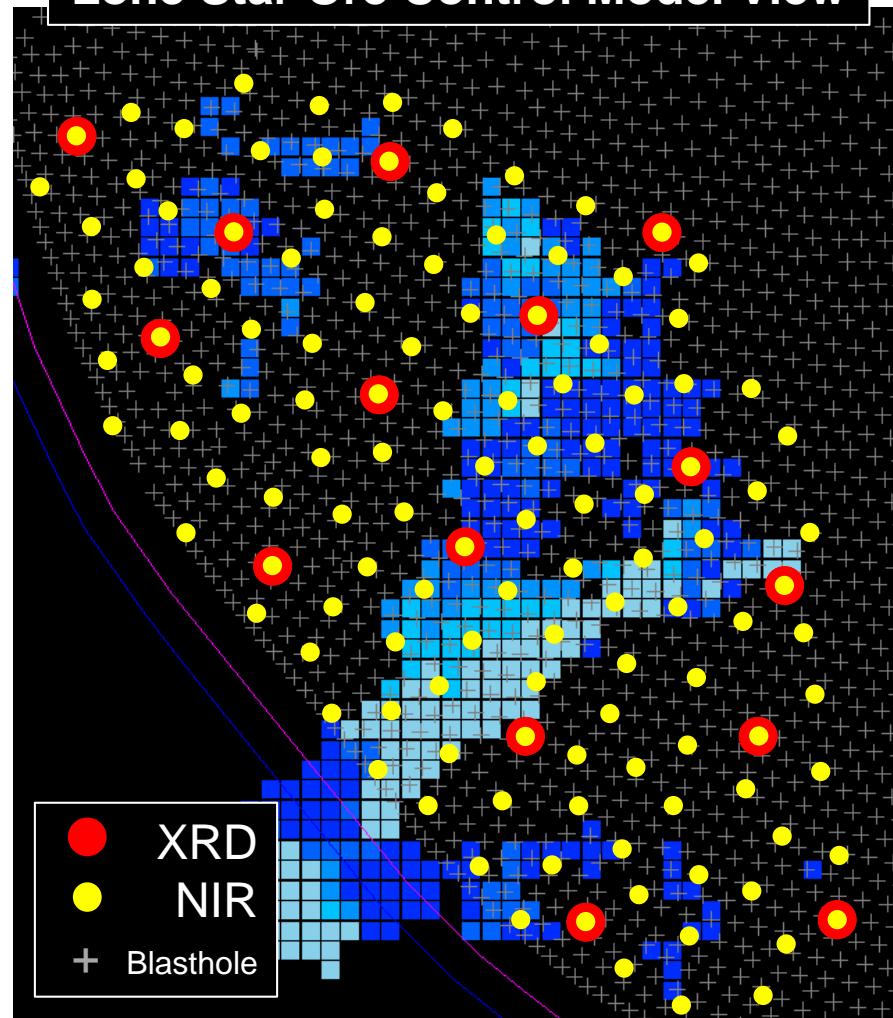
- **50%** of blastholes sampled for NIR, **20%** of blastholes sampled for XRD

- NIR data used to calculate **total clay**

Total Clay = Kaolinite (%) + Swelling Clay (%)

- Total clay used to determine if **blending** is necessary

Lone Star Ore Control Model View



XRD is ~8x the cost of NIR per sample

Safford: Direct Application of NIR – Crush/Leach Circuit

Effective Clay Blending:

- Prevents **clogging** of grizzly feeder in crusher
- Allows for more **effective agglomeration** stage



Safford: Direct Application of NIR - Leachpad

Effective Clay Blending:

- Allows for **effective percolation** of sulfuric acid down through lifts
- Prevents **geotechnical instabilities** in leach pad construction



Questions?