



nimbus
INFORMATICS

<http://nimbusinformatics.com>

<http://lifetechnologies.com/>

The import of sequence data on a private cluster install of SeqWare is straightforward. However, maintaining IRB and HIPAA compliance on public clouds such as EC2 requires careful security procedures. SeqWare provides tools to facilitate the encryption of data into and out of the cloud. This includes encryption of sequence data sent by users and results sent back, encrypted network connections and filesystems on compute nodes, and following Linux security best practices.

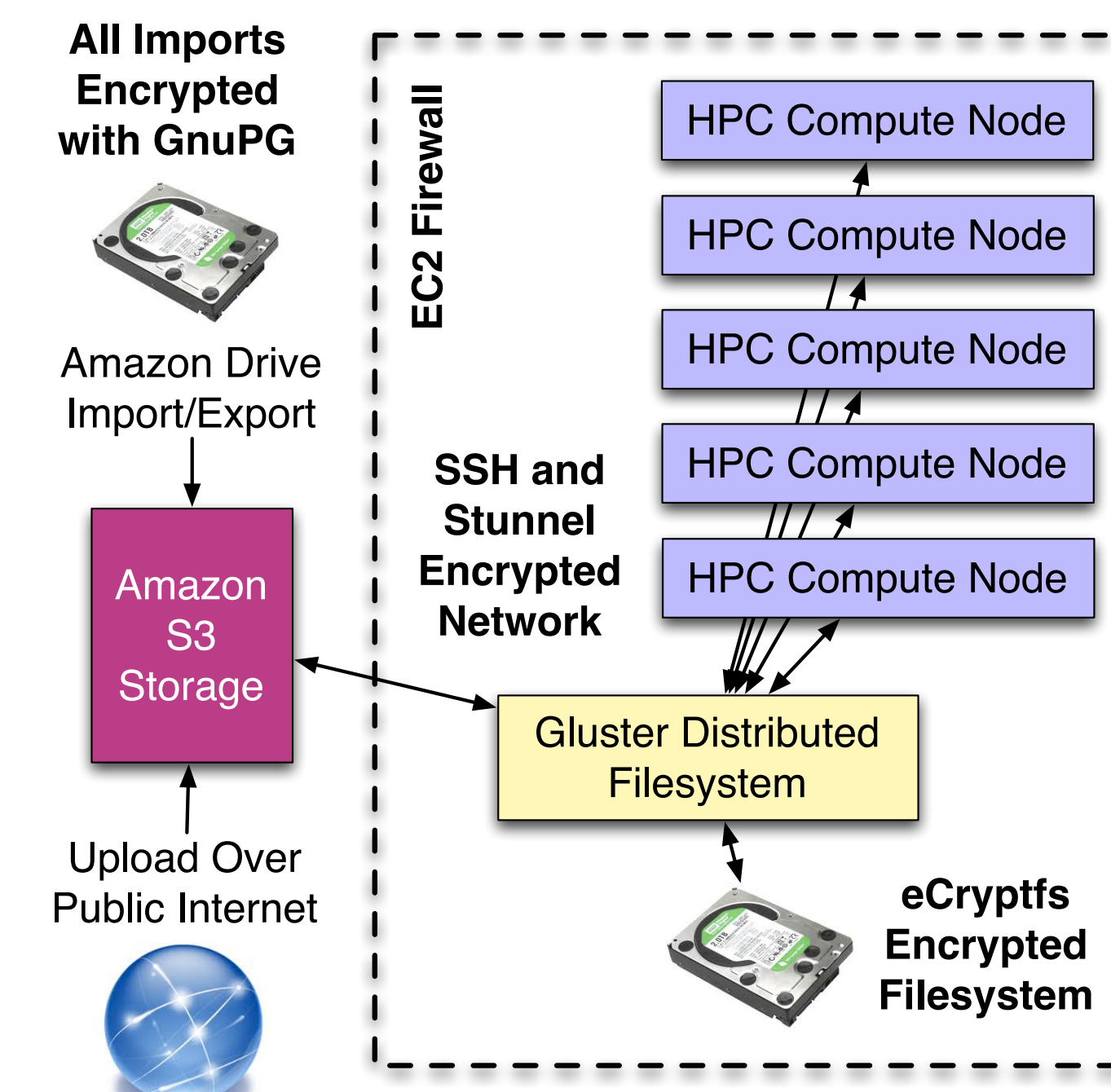


Figure 3: Security in the Amazon Cloud

Genome	Step	Current Timing	Current Cost	Projected Timing	Projected Costs	Item	Period	Cost
	Data Import (drive import)	~2-4 days	\$100	~2-4 days	\$100	Recurring Storage Costs	per month	\$5
	Temp Data Storage	~1 week	\$300	~3 days	\$100	Recurring Archive Costs	per month	\$45
	HPC Computation Time	5 days, 8 nodes	\$1,500	3 days, 8 nodes	\$900	Recurring Hosting Costs	per month	\$15
	Total	~8 days	\$1,900	~6 days	\$1,100	Total	per month	\$65
Exome	Step	Current Timing	Current Cost	Projected Timing	Projected Costs	Item	Period	Cost
	Data Import (upload)	~1 days	\$10	~1 days	\$10	Recurring Storage Costs	per month	\$5
	Temp Data Storage	~3 days	\$30	~1 days	\$10	Recurring Archive Costs	per month	\$5
	HPC Computation Time	2 days, 2 nodes	\$150	1 days, 2 nodes	\$75	Recurring Hosting Costs	per month	\$15
	Total	~2 days	\$190	~1 day	\$95	Total	per month	\$25

Workflow Automation

[illegible]

Figure 4: The Lucier genome browser and query engine are available at <http://lucier.nimbusinformatics.com>

Massively parallel sequencing technologies have opened the door to a future where thousands of genomes will be sequenced on a regular basis. A considerable challenge exists for how to analyze and query the massive amount of data produced. Here we have presented SeqWare, a flexible open source framework designed to run both on a local environment and the cloud, as a potential solution. The project is fully open source and available from <http://seqware.sf.net> for deployment locally or on EC2. Nimbus Informatics LLC provides a commercially-supported, easy-to-use web-based version of SeqWare backed by Amazon's EC2 cloud (currently in private beta, please see <http://nimbusinformatics.com>). SeqWare development is supported by Nimbus Informatics LLC, Life Technologies Inc., users including UNC, UCLA, and the Cancer Genome Atlas, and the open source community.

2 Nimbus Informatics, LLC sponsors the open source SeqWare project and provides commercial support, data hosting, and push-button utilization of SeqWare through an easy-to-use web interface. This service is currently in private beta, see <http://nimbusinformatics.com> for more information.