

Document, Store, Retrieve and Share your Science

- Create, manage and share experiments and notebooks with your colleagues, with access to the new scientific collaboration platform, PerkinElmer Signals™ Notebook Individual Edition
- Draw reaction schemes using a web enabled ChemDraw sketcher, the stoichiometry is calculated for you
- Add data from any source, of any type, from any browser
- Free text and chemical search and query help you find your data, whenever, wherever
- A cloud based solution means no installation, frequent updates, seamless upgrades, no migrations – no overhead

Analyze and Explore Your Reactions

- Identify similar reactions from the public literature with integration with SciFinder (Additional SciFinder license required)
- Perform enhanced retrosynthesis analyses on molecules to identify the reactions that are required to create a bond or set of bonds emanating from a single atom
- Easily manage the numbering of your molecules with dynamic Reaction Auto-numbering
- Automatically calculate, track and update stoichiometry data for chemical reactions

Predict and Calculate Efficiently and Accurately – Every Time

- Predict and calculate values for physicochemical properties important for bioavailability and stability, such as mp, bp, cLogP, pKa, LogD, and LogS
- Predict ¹H and ¹³C 1D NMR spectra, including varying solvent and frequency, update with your own experimental data to improve the predictivity
- Explore the 3D shape and properties of compounds with the Chem3D® molecular modeling and protein visualization tool
- Add chemical intelligence to Microsoft® Excel® spreadsheets. Build and manipulate chemical structures, compute chemical properties and use structure and substructure searches to locate and group compounds
- Extend the capabilities of ChemDraw by developing your own JavaScript-based add-ins.

ChemDraw Everywhere

- Store, retrieve and edit your ChemDraw documents on the go, using ChemDraw Cloud
- Web based ChemDraw sketcher allows access from any device
- All documents are stored, encrypted, and backed-up securely on ChemDraw Cloud servers
- Embed ChemDraw into your custom web applications with ChemDraw JavaScript (with ChemOffice Professional Site License)

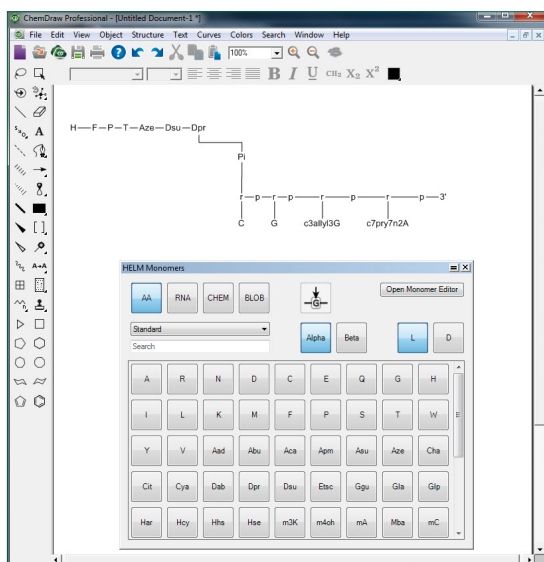


Figure 3: Create novel, complex biopolymers with the new HELM editor.