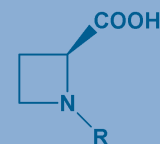
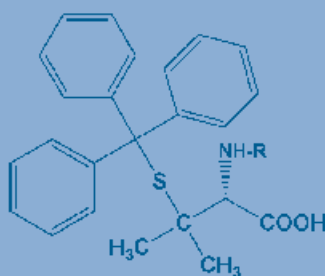
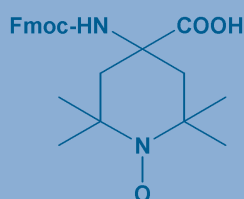
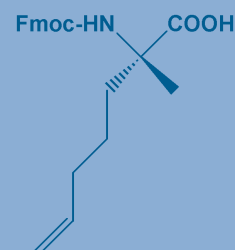
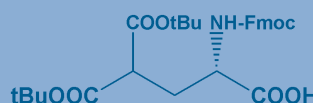
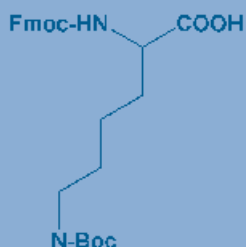
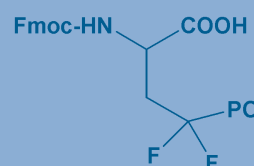
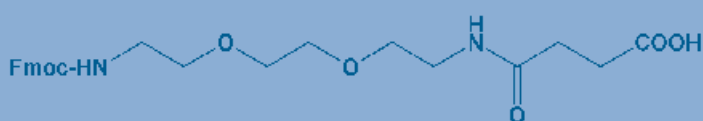


Large collections

amino acids

building blocks / unusual amino acids



Certified to ISO 9001:2008

