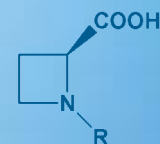
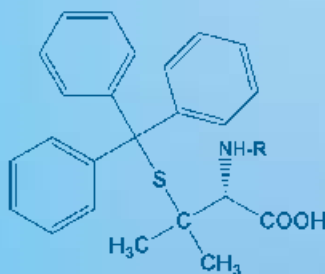
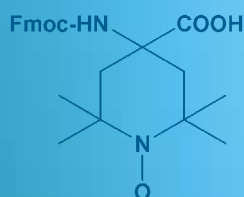
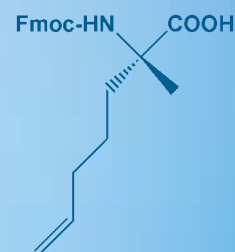
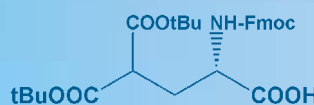
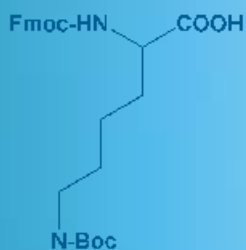
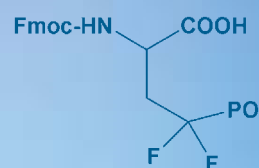
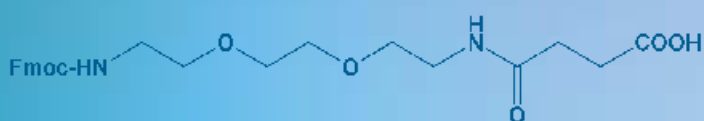


L a r g e c o l l e c t i o n s

# amino acids

building blocks / unusual amino acids



ISO 9001:2000 Company

