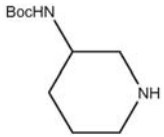
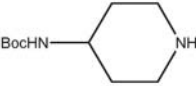
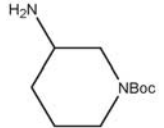
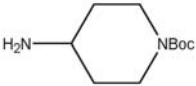

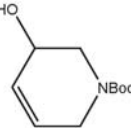
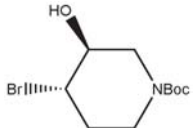


# New Piperidine Building Blocks

Product	Structure	Size	Catalog #	US\$
3-tert-Butoxycarbonyl-aminopiperidine		1 g 5 g	60010-1 60010-5	\$150 \$600
4-tert-Butoxycarbonyl-aminopiperidine		5 g 25 g	60011-5 60011-25	\$120 \$565
N-1-Boc-3-aminopiperidine		1 g 5 g	60012-1 60012-5	\$120 \$480
N-1-Boc-4-aminopiperidine		1 g 5 g	60013-1 60013-5	\$105 \$420
N-Boc-1,2,3,6-tetrahydropyridine		1 g 5 g	60014-1 60014-5	\$80 \$320
N-Boc-3-hydroxy-1,2,3,6-tetrahydropyridine		0.25 g 1 g 5 g	60015-025 60015-1 60015-5	\$75 \$200 \$800
N-Boc-trans-4-bromo-3-hydroxypiperidine		1 g 5 g	60016-1 60016-5	\$125 \$500