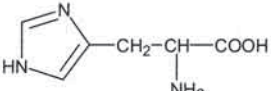
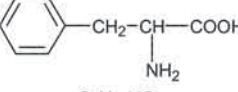
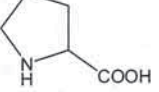
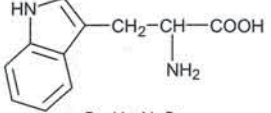
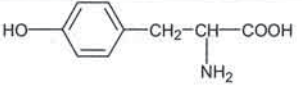


参考資料 6

アミノ酸の構造・略号 (3文字, 1文字)

$\begin{array}{c} \text{H}_3\text{C}-\text{CH}-\text{COOH} \\ \\ \text{NH}_2 \end{array}$ <p>C₃H₇NO₂ MW = 89.09 (71.07)</p> <p>Alanine (アラニン) Ala (A)</p>	$\begin{array}{c} \text{H}_2\text{N}-\text{C}-\text{NH}-(\text{CH}_2)_3-\text{CH}-\text{COOH} \\ \qquad \qquad \\ \text{NH} \qquad \qquad \text{NH}_2 \end{array}$ <p>C₆H₁₄N₄O₂ MW = 174.20 (156.18)</p> <p>Arginine (アルギニン) Arg(R)</p>	$\begin{array}{c} \text{H}_2\text{N}-\text{C}-\text{CH}_2-\text{CH}-\text{COOH} \\ \qquad \qquad \\ \text{O} \qquad \qquad \text{NH}_2 \end{array}$ <p>C₄H₉N₂O₃ MW = 132.12 (114.10)</p> <p>Asparagine (アスパラギン) Asn(N)</p>
$\begin{array}{c} \text{HOOC}-\text{CH}_2-\text{CH}-\text{COOH} \\ \\ \text{NH}_2 \end{array}$ <p>C₄H₇NO₄ MW = 133.10 (115.08)</p> <p>Asparatic acid (アスパラギン酸) Asp(D)</p>	$\begin{array}{c} \text{HS}-\text{CH}_2-\text{CH}-\text{COOH} \\ \\ \text{NH}_2 \end{array}$ <p>C₃H₇NO₂S MW = 121.16 (103.14)</p> <p>Cysteine (システイン) Cys(C)</p>	$\begin{array}{c} \text{H}_2\text{N}-\text{C}-\text{CH}_2-\text{CH}-\text{COOH} \\ \qquad \qquad \\ \text{O} \qquad \qquad \text{NH}_2 \end{array}$ <p>C₅H₁₀N₂O₃ MW = 146.14 (128.13)</p> <p>Glutamine (グルタミン) Gln(Q)</p>
$\begin{array}{c} \text{HOOC}-(\text{CH}_2)_2-\text{CH}-\text{COOH} \\ \\ \text{NH}_2 \end{array}$ <p>C₅H₉NO₄ MW = 147.13 (129.11)</p> <p>Glutamic acid (グルタミン酸) Glu(E)</p>	$\begin{array}{c} \text{H}-\text{CH}-\text{COOH} \\ \\ \text{NH}_2 \end{array}$ <p>C₂H₅NO₂ MW = 75.07 (57.05)</p> <p>Glycine (グリシン) Gly (G)</p>	 <p>C₆H₉N₃O₂ MW = 155.15 137.14</p> <p>Histidine (ヒスチジン) His(H)</p>
$\begin{array}{c} \text{H}_3\text{C}-\text{CH}_2 \\ \qquad \\ \text{CH}-\text{CH}-\text{COOH} \\ \qquad \\ \text{H}_3\text{C} \qquad \text{NH}_2 \end{array}$ <p>C₆H₁₃NO₂ MW = 131.17 (113.15)</p> <p>Isoleucine (イソロイシン) Ile (I)</p>	$\begin{array}{c} \text{H}_3\text{C} \\ \qquad \\ \text{CH}-\text{CH}_2-\text{CH}-\text{COOH} \\ \qquad \\ \text{H}_3\text{C} \qquad \text{NH}_2 \end{array}$ <p>C₆H₁₃NO₂ MW = 131.17 (113.15)</p> <p>Leucine (ロイシン) Leu (L)</p>	$\begin{array}{c} \text{H}_2\text{N}-(\text{CH}_2)_4-\text{CH}-\text{COOH} \\ \\ \text{NH}_2 \end{array}$ <p>C₆H₁₄N₂O₂ MW = 146.19 (128.17)</p> <p>Lysine (リシン) Lys(K)</p>
$\begin{array}{c} \text{H}_3\text{C}-\text{S}-(\text{CH}_2)_2-\text{CH}-\text{COOH} \\ \\ \text{NH}_2 \end{array}$ <p>C₆H₁₁NO₂S MW = 149.21 (131.19)</p> <p>Methionine (メチオニン) Met(M)</p>	 <p>C₉H₉NO₂ MW = 165.19 (147.17)</p> <p>Phenylalanine (フェニルアラニン) Phe(F)</p>	 <p>C₅H₉NO₂ MW = 115.13 (97.11)</p> <p>Proline (プロリン) Pro(P)</p>
$\begin{array}{c} \text{HO}-\text{CH}_2-\text{CH}-\text{COOH} \\ \\ \text{NH}_2 \end{array}$ <p>C₃H₇NO₃ MW = 105.09 (87.07)</p> <p>Serine (セリン) Ser (S)</p>	$\begin{array}{c} \text{HO} \\ \\ \text{CH}-\text{CH}-\text{COOH} \\ \qquad \\ \text{H}_3\text{C} \qquad \text{NH}_2 \end{array}$ <p>C₄H₉NO₃ MW = 119.12 (101.10)</p> <p>Threonine (トレオニン) Thr (T)</p>	 <p>C₁₁H₁₂N₂O₂ MW = 204.23 (186.20)</p> <p>Tryptophan (トリプトファン) Tyr(W)</p>
 <p>C₉H₉NO₃ MW = 181.19 (163.17)</p> <p>Tyrosine (チロシン) Tyr(Y)</p>	$\begin{array}{c} \text{H}_3\text{C} \\ \qquad \\ \text{CH}-\text{CH}-\text{COOH} \\ \qquad \\ \text{H}_3\text{C} \qquad \text{NH}_2 \end{array}$ <p>C₅H₁₁NO₂ MW = 117.15 (99.13)</p> <p>Valine (バリン) Val (V)</p>	<p><表記例></p> <p>$\begin{array}{c} \text{H}_3\text{C}-\text{CH}-\text{COOH} \\ \\ \text{NH}_2 \end{array}$ ← 構造式</p> <p>C₃H₇NO₂ MW = 89.09 (71.07) ← 分子量 分子量 (残基量)</p> <p>Alanine (アラニン) Ala (A) ← 名称 略号</p>

※残基量：アミノ酸分子量より水の分子量を差し引いたもの