

Benzamil . HCl

Na⁺ channel blocker

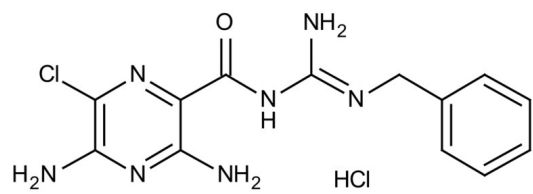
Derivative of amiloride (Prod. No. BML-CA200). Potent sodium channel blocker.

Citations: 5

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

| | |
|--------------------------|--|
| Use/Stability | As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at room temperature. Stock solutions are stable for up to 3 months at -20°C. |
| Handling | Protect from light. |
| Long Term Storage | Ambient |
| Shipping | Ambient Temperature |

Regulatory Status RUO - Research Use Only

Product Details

| | |
|-------------------------|--|
| Alternative Name | Benzylamiloride . HCl, N-(Benzylamidino)-3,5-diamino-6-chloropyrazinecarboxamide . HCl |
| Appearance | Yellow solid. |
| CAS | 161804-20-2 |
| Couple Target | Sodium channel |
| Couple Type | Blocker |
| Formula | $C_{13}H_{14}ClN_7O$. HCl |
| MW | 356.2 |
| Purity | ≥98% (HPLC) |
| Solubility | Soluble in water. |

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)